



# 2024 PRODUCT CATALOG





**SignalChem is present & active on  
these social media channels**



<https://www.linkedin.com/company/signalchem>



<https://www.facebook.com/SignalChem/>

## **Sign up for our Newsletters!**

We send informative, engaging, and educational newsletters



For more information, please write us on:  
[info@signalchem.com](mailto:info@signalchem.com)

# Contents

## Enzymes

### Kinases

Protein Kinases

Lipid Kinases

Carbohydrate Kinases

Mutant Kinases

### Phosphatases

### Epigenetic Enzymes

Histone Deacetylases

Acetyl & Methyltransferases

Deacetylase & Demethylase Proteins

Bromodomains

### Ubiquitins

Ubiquitin Enzymes - E1

Ubiquitin Enzymes - E2

Ubiquitin Enzymes - E3

Ubiquitin Enzymes - DUBs

Ubiquitin-like Proteins

### Arginine Deiminase & Proteins

### G Proteins

GTPases

Other G-Proteins

GAPs/GEFs

### Proteases

### Phosphodiesterases

### Other Enzymes

Oxidoreductases

Hydrolases

Transferases

## Antibodies

Isoform Specific Antibodies

Tag Antibodies

Primary Antibodies

Secondary Antibodies

Modification-Specific Antibodies

Phospho-Specific Antibodies

## Proteins

Adaptor Proteins

Apoptosis

Cell-Cycle Proteins

Cell Stress and Chaperone Proteins

Fluorescent and Color Proteins

Microtubule and Actin-Associated Proteins

Regulatory Proteins

Transcription Proteins

Tau Proteins

Growth Factors

Affinity Proteins and Resins

## Compounds

Epigenetic Enzyme Activators

Kinase Inhibitors

Phosphatase Inhibitors

Epigenetic Enzyme Inhibitors

Kinase Activators

## Reagents

Assay Buffer and Cofactors

Assay Substrates

Lipid Substrates

Oligo Substrates

Carbohydrate Substrates

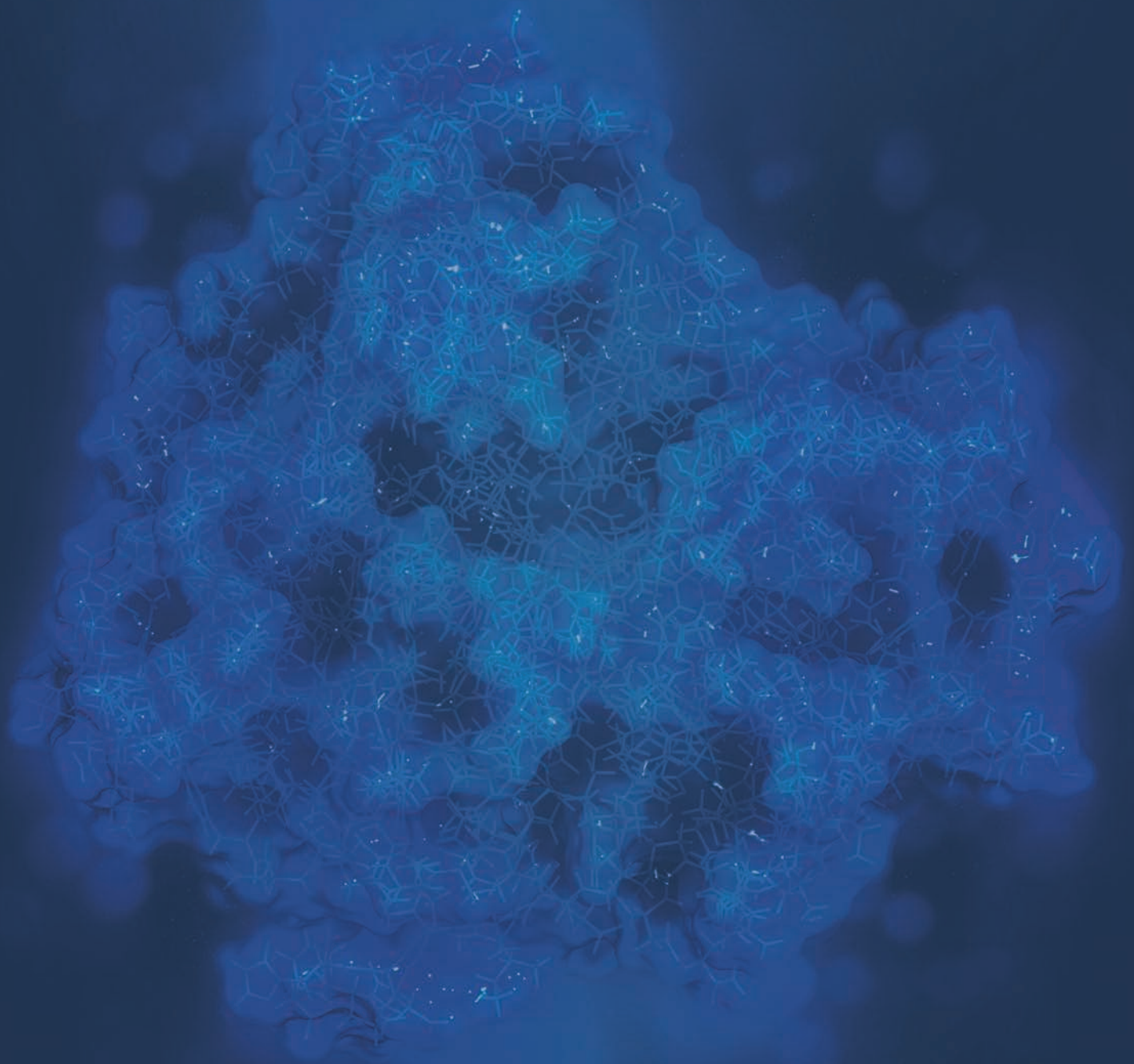
Peptide Substrates

Protein Substrates

## PAD Kits

Detection kit

Inhibitor Screening Kit



World's Largest  
**Kinase**  
Catalog

# Enzymes

## Kinases

### Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
AAK1, Active	A01-11G	Human	GST	Sf9 Cells	1-465
AATK Protein	A02-35G	Mouse	GST	Sf9 Cells	8-411
ABL1, Active	A03-11H	Mouse	HIS	Sf9 Cells	27-end
ABL1, Active	A03-18H	Human	HIS	Sf9 Cells	27-end
ABL2, Active	A04-11H	Human	HIS	Sf9 Cells	38-end
ACK, Active	A05-11G	Human	GST	Sf9 Cells	110-476
ACKA Protein	A120-30H	Bacteria	HIS	E.coli	Full Length
ACKA Protein	A120-30G	Bacteria	GST	E.coli	Full Length
ACTR2 Protein	A39-30G	Human	GST	Sf9 Cells	Full Length
ACVR2A, Active	A35-11H	Human	HIS	Sf9 Cells	161-end
ACVR2B, Active	A36-11G	Human	GST	Sf9 Cells	182-488
ADCK1 Protein	A10-11G	Human	GST	Sf9 Cells	18-end
ADCK3 Protein	A37-11G	Human	GST	Sf9 Cells	156-end
ADCK5 Protein	A13-11G	Human	GST	Sf9 Cells	68-end
AKT1, Active	A16-10G	Human	GST	Sf9 Cells	Full Length
AKT1, Active	A16-10H	Human	HIS	Sf9 Cells	Full-length
AKT1, Unactive	A16-14G	Human	GST	Sf9 Cells	Full Length
AKT2, Active	A17-10H	Human	HIS	Sf9 Cells	Full Length
AKT2, Active	A17-10G	Human	GST	Sf9 Cells	Full Length
AKT2, Unactive	A17-15H	Human	HIS	Sf9 Cells	142-end
AKT2, Unactive	A17-14G	Human	GST	Sf9 Cells	Full Length
AKT3, Active	A18-10G	Human	GST	Sf9 Cells	Full Length
AKT3, Unactive	A18-14G	Human	GST	Sf9 Cells	Full Length
ALK, Active	A19-11G	Human	GST	Sf9 Cells	1060-end
ALK1, Active	A09-11G	Human	GST	Sf9 Cells	144-end
ALK2, Active	A06-11G	Human	GST	Sf9 Cells	147-end
ALK2, Active	A06M-11G	Mouse	GST	Sf9 Cells	147-end
ALK3 (BMPR1A), Active	B04-11G	Human	GST	Sf9 Cells	187-end
ALK4, Active	A07-11G	Human	GST	Sf9 Cells	150-end
ALK6 (BMPR1B), Active	B05-11G	Human	GST	Sf9 Cells	149-end
ALK7 Protein	A08-35G	Human	GST	Sf9 Cells	135-end
AMHR2 Protein	A38-35G	Human	GST	Sf9 Cells	171-end
AMPK (A1/B1/G1), Active	P47-10H	Human	HIS	Sf9 Cells	Full Length
AMPK (A1/B1/G1), Active	P47-110GH	Human	GST/HIS	Sf9 Cells	Full length
AMPK (A1/B1/G2), Active	P55-10H	Human	HIS	Sf9 Cells	Full Length
AMPK (A1/B1/G2), Active	P55-10G	Human	GST/HIS	Sf9 Cells	Full Length
AMPK (A1/B1/G3), Active	P56-10H	Human	HIS	Sf9 Cells	Full Length
AMPK (A1/B1/G3), Active	P56-10G	Human	GST/HIS	Sf9 Cells	Full Length
AMPK (A1/B2/G1), Active	P50-10H	Human	HIS	Sf9 Cells	Full Length
AMPK (A1/B2/G1), Active	P50-10G	Human	GST/HIS	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
AMPK (A1/B2/G2), Active	P45-10G	Human	GST/HIS	Sf9 Cells	Full Length
AMPK (A1/B2/G3), Active	P83-10G	Human	GST/HIS	Sf9 Cells	Full Length
AMPK (A2/B1/G1), Active	P48-10H	Human	HIS	Sf9 Cells	Full Length
AMPK (A2/B1/G1), Unactive	P48-14H	Human	HIS	Sf9 Cells	Full Length
AMPK (A2/B1/G2), Active	P84-10GH	Human	GST/HIS	Sf9 Cells	Full Length
AMPK (A2/B1/G3), Active	P85-10GH	Human	GST/HIS	Sf9 Cells	Full Length
AMPK (A2/B2/G1), Active	P49-10H	Human	HIS	Sf9 Cells	Full Length
AMPK (A2/B2/G1), Active	P49-10GH	Human	GST/HIS	Sf9 Cells	Full Length
AMPK (A2/B2/G1), Active	P49M-10H	Mouse	HIS	Sf9 Cells	Full-Length
AMPK (A2/B2/G2), Active	P54-10H	Human	HIS	Sf9 Cells	Full Length
AMPK (A2/B2/G2), Active	P54-10GH	Human	GST/HIS	Sf9 Cells	Full Length
AMPK (A2/B2/G3), Active	P46-10GH	Human	GST/HIS	Sf9 Cells	Full Length
AMPK (A2/B2/G3), Unactive	P46-14GH	Human	GST/HIS	Sf9 Cells	Full Length
AMPK(A1/B2/G1), Active	P50M-10H	Mouse	HIS	Sf9 Cells	Full Length
ANKK1 (SGK288), Active	A20-10G	Human	GST	Sf9 Cells	Full Length
ARAF Protein	A24-35G	Human	GST	Sf9 Cells	209-end
ASK1, Active	M13-11G	Human	GST	Sf9 Cells	649-946
ASK1, Active	M13-10G	Human	GST	Sf9 Cells	Full Length
ASK3 Protein	M73-31G	Human	GST	Sf9 Cells	582-1169
ASK3, Active	M73-10G	Human	GST	Sf9 Cells	Full-Length
ATM Protein	A26-35G	Human	GST	Sf9 Cells	2667-end
ATR Protein	A27-35G	Human	GST	Sf9 Cells	2245-end
AURORA A, Active	A28-10G	Mouse	GST	Sf9 Cells	Full Length
AURORA A, Active	A28-18G	Human	GST	Sf9 Cells	Full Length
AURORA A, Active	A28-18H	Human	HIS	Sf9 Cells	Full Length
AURORA A, Unactive	A28-14G	Human	GST	Sf9 Cells	Full Length
AURORA B, Active	A31-10G	Human	GST	Sf9 Cells	Full Length
AURORA B, Active	A31-10H	Human	HIS	Sf9 Cells	Full Length
AURORA B, Unactive	A31-14G	Human	GST	Sf9 Cells	Full Length
AURORA C, Active	A33-10G	Human	GST	Sf9 Cells	Full Length
AURORA C, Unactive	A33-14G	Human	GST	Sf9 Cells	Full Length
AXL, Active	A34-11H	Human	HIS	Sf9 Cells	473-end
AXL, Active	A34-11G	Human	GST	Sf9 Cells	473-end
BCKDK Protein	B15-31G	Human	GST	Sf9 Cells	30-end
BCR Protein	B01-35G	Human	GST	Sf9 Cells	1-695
BLK, Active	B02-10G	Human	GST	Sf9 Cells	Full Length
BMP2K, Active	B03B-11G	Mouse	GST tag	Sf9 Cells	1-422
BMPR2, Active	B06-11H	Human	HIS	Sf9 Cells	174-end
BMX, Active	B07-10G	Human	GST	Sf9 Cells	Full Length
BMX, Active	B07-10CG	Rat	GST	Sf9 Cells	Full Length
BRAF, Active	B08-11G	Human	GST	Sf9 Cells	416-766
BRAF, Active	B08-11BG	Human	GST	Sf9 Cells	381-end or exon 9-18
BRK, Active	P94-10G	Human	GST	Sf9 Cells	Full Length
BRSK1, Active	B13-10G	Mouse	GST	Sf9 Cells	Full Length
BRSK2, Active	B14-10G	Human	GST	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
BTK, Active	B10-10H	Human	HIS	Sf9 Cells	Full Length
BTK, Active	B10-10CG	Rat	GST	Sf9 Cells	Full Length
BUB1 Protein	B11-35G	Human	GST	Sf9 Cells	756-end
BUB1B, Active	B12-11G	Human	GST	Sf9 Cells	613-end
c-KIT, Active	K06-11BG	Human	GST	Sf9 Cells	544-end
c-MER, Active	M51-11G	Human	GST	Sf9 Cells	578-872
c-MER, Active	M51-112G	Human	GST	Sf9 Cells	528-end
c-MER, Active	M51-172G	Mouse	GST	Sf9 Cells	528-end
CAMK1 beta, Active	C08-10G	Mouse	GST	Sf9 Cells	Full Length
CAMK1 delta, Active	C09-10G	Human	GST	Sf9 Cells	Full Length
CAMK1 gamma, Active	C10-11G	Human	GST	Sf9 Cells	1-330
CAMK1 gamma, Active	C10-10BG	Human	GST	Sf9 Cells	Full Length
CAMK1, Active	C07-10G	Mouse	GST	E.coli	Full Length
CAMK2 alpha, Active	C11-10G	Human	GST	Sf9 Cells	Full Length
CAMK2 beta, Active	C12-10H	Human	HIS	Sf9 Cells	Full Length
CAMK2 beta, Active	C12-10G	Human	GST	Sf9 Cells	Full Length
CAMK2 delta, Active	C13-10G	Human	GST	Sf9 Cells	Full Length
CAMK2 gamma, Active	C14-10G	Human	GST	Sf9 Cells	c-terminal truncation
CAMK2 gamma, Active	C14-10BG	Human	GST	Sf9 Cells	Full Length
CAMK4, Active	C15-10G	Human	GST	Sf9 Cells	Full Length
CAMKK1, Active	C17-10G	Human	GST	Sf9 Cells	Full Length
CAMKK1, Active	C17-18G	Rat	GST	Sf9 Cells	Full Length
CAMKK2, Active	C18-10G	Human	GST	Sf9 Cells	Full Length
CASK, Active	C19-11G	Human	GST	Sf9 Cells	1-570
CDC42BPG (DMPK2), Active	D17-11G	Human	GST	Sf9 Cells	1-475
CDC7/DBF4, Active	C26-10G	Human	GST	Sf9 Cells	Full Length
CDK1, Unactive	C22-14G	Human	GST	Sf9 Cells	Full Length
CDK1/CyclinA1, Active	C22-18BG	Human	GST	Sf9 Cells	Full Length
CDK1/CyclinA2, Active	C22-18G	Human	GST	Sf9 Cells	Full Length
CDK1/CyclinB1, Active	C22-10G	Human	GST	Sf9 Cells	Full Length
CDK10/CyclinM Protein	C41-34G	Human	GST	Sf9 Cells	Full length
CDK11, Unactive	C42-14G	Human	GST	Sf9 Cells	Full Length
CDK11/CyclinC Protein	C42-34G	Human	GST	Sf9 Cells	Full Length
CDK11A (CDC2L2) Protein	C24-31G	Human	GST	Sf9 Cells	361-end
CDK11A (CDC2L2)/CyclinD3 Protein	C24-35CG	Human	GST	Sf9 Cells	361-end (CDK11A) and Full Length (CyclinD3)
CDK11B (CDC2L1) Protein	C23-30G	Human	GST	Sf9 Cells	Full Length
CDK12 (CRKRS) Protein	C62-35G	Human	GST	Sf9 Cells	658-end
CDK12/CyclinK, Active	C62-11BG	Human	GST	Sf9 Cells	Full Length
CDK13 (CDC2L) Protein	C25-35G	Human	GST	Sf9 Cells	516-end
CDK13/CyclinK, Active	C25-11BG	Human	GST	Sf9 Cells	Full length
CDK2, Unactive	C29-14BG	Human	GST	Sf9 Cells	Full Length
CDK2/CyclinA1, Active	C29-10BG	Human	GST	Sf9 Cells	Full Length
CDK2/CyclinA2, Active	C29-10G	Human	GST	Sf9 Cells	Full Length
CDK2/CyclinE1, Active	C29-18G	Human	GST	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
CDK2/CyclinE1, Unactive	C29-14G	Human	GST	Sf9 Cells	Full Length
CDK2/CyclinE2, Active	C29-18BG	Human	GST	Sf9 Cells	Full Length
CDK2/CyclinO, Active	C29-19G	Human	GST	Sf9 Cells	Full Length
CDK20 (CCRK) Protein	C21-30G	Mouse	GST	Sf9 Cells	Full Length
CDK3, Unactive	C30-14G	Human	GST	Sf9 Cells	Full Length
CDK3/CyclinE1, Active	C30-10G	Human	GST	Sf9 Cells	Full Length
CDK3/CyclinE2, Active	C30-10BG	Human	GST	Sf9 Cells	Full Length
CDK4, Unactive	C31-14G	Human	GST	E.coli	Full Length
CDK4, Unactive	C31-14BG	Human	GST	Sf9 Cells	Full Length
CDK4/CyclinD1, Active	C31-10G	Human	GST	Sf9 Cells	Full Length
CDK4/CyclinD3, Active	C31-18G	Human	GST	Sf9 Cells	Full Length
CDK5, Unactive	C33-14G	Human	GST	Sf9 Cells	Full Length
CDK5/p25, Active	C33-10G	Human	GST	Sf9 Cells	Full Length
CDK5/p35, Active	C33-10BG	Human	GST	Sf9 Cells	Full Length
CDK6, Unactive	C35-15H	Human	HIS	Sf9 Cells	Full Length
CDK6/CyclinD1, Active	C35-18H	Human	GST/HIS	Sf9 Cells	Full Length
CDK6/CyclinD2 Protein	C35-34G	Human	GST	Sf9 Cells	Full Length
CDK6/CyclinD3, Active	C35-10H	Human	HIS	Sf9 Cells	Full Length
CDK6/CyclinD3, Unactive	C35-14H	Human	HIS	Sf9 Cells	Full Length
CDK7, Unactive	C36-14G	Human	GST	Sf9 Cells	Full Length
CDK7/CyclinH1, Active	C36-102H	Human	HIS	Sf9 Cells	Full Length
CDK7/CyclinH1/MNAT1, Active	C36-10H	Human	HIS	Sf9 Cells	Full Length
CDK8, Unactive	C38-14G	Human	GST	Sf9 Cells	Full Length
CDK8/CyclinC Protein	C38-34G	Human	GST	Sf9 Cells	Full Length
CDK9, Unactive	C40-14G	Human	GST	Sf9 Cells	Full Length
CDK9/CyclinK, Active	C40-10G	Human	GST	Sf9 Cells	Full Length
CDK9/CyclinT1, Active	C40-18CG	Human	GST	Sf9 Cells	Full Length
CDK9/CyclinT2, Active	C40-18BG	Human	GST	Sf9 Cells	Full Length
CDKL1 Protein	C43-30G	Human	GST	Sf9 Cells	Full Length
CDKL2 Protein	C44-30G	Human	GST	Sf9 Cells	Full Length
CDKL3 Protein	C45-30G	Human	GST	Sf9 Cells	Full Length
CDKL4 Protein	C69-30G	Human	GST	Sf9 Cells	Full Length
CHK1, Active	C47-10G	Human	GST	Sf9 Cells	Full Length
CHK1, Active	C47-10H	Human	HIS	Sf9 Cells	Full Length
CHK2, Active	C48-10G	Human	GST	Sf9 Cells	Full Length
CHK2, Active	C48-10H	Human	HIS	Sf9 Cells	Full length
CK1 alpha 1, Active	C64-10G	Human	GST	Sf9 Cells	Full Length
CK1 alpha 1L, Active	C75-10G	Human	GST	Sf9 Cells	Full Length
CK1 delta, Active	C65-10G	Human	GST	Sf9 Cells	Full Length
CK1 epsilon, Active	C66-10G	Human	GST	Sf9 Cells	Full Length
CK1 gamma 1, Active	C68-11G	Human	GST	Sf9 Cells	21-end
CK1 gamma 2, Active	C68-10BG	Human	GST	Sf9 Cells	Full Length
CK1 gamma 3, Active	C68-10CG	Human	GST	Sf9 Cells	Full Length
CK2 alpha 1, Active	C70-10G	Human	GST	Sf9 Cells	Full Length
CK2 alpha 2, Active	C71-10G	Human	GST	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
CLIK1L Protein	C76-34G	Human	GST	Sf9 Cells	Full-Length
CLK1, Active	C57-11G	Human	GST	Sf9 Cells	129-end
CLK2, Active	C58-11G	Human	GST	Sf9 Cells	137-end
CLK3, Active	C60-10G	Human	GST	Sf9 Cells	Full Length
CLK4, Active	C61-11G	Human	GST	Sf9 Cells	70-end
COT, Active	M16-11G	Human	GST	Sf9 Cells	30-397
COT, Unactive	M16-15G	Human	GST	Sf9 Cells	30-397
CSK, Active	C63-10G	Human	GST	Sf9 Cells	Full Length
CSK, Active	C63-10U	Human	TAG-Free	Sf9 Cells	Full Length
DAPK1, Active	D01-11G	Human	GST	Sf9 Cells	1-363
DAPK2, Active	D02-10G	Human	GST	Sf9 Cells	Full Length
DAPK3, Active	D03-10G	Human	GST	Sf9 Cells	Full Length
DCAMKL1, Active	D14-10G	Human	GST	Sf9 Cells	Full Length
DCAMKL1, Active	D14-11G	Human	GST	Sf9 Cells	266-end
DCAMKL1, Active	D14-10H	Human	HIS	Sf9 Cells	Full Length
DCAMKL2, Active	D15-11G	Human	GST	Sf9 Cells	1-690
DDR1, Active	D05-11G	Human	GST	Sf9 Cells	444-end
DDR2, Active	D06-11BG	Human	GST	Sf9 Cells	467-end
DDR2, Active	D06M-11BG	Mouse	GST	Sf9 Cells	467-end
DLK, Active	M20-11G	Human	GST	Sf9 Cells	1-520
DMPK, Active	D07-10G	Human	GST	Sf9 Cells	Full Length
DNAPK Protein	D08-35G	Human	GST	Sf9 Cells	3746-end
DRAK1 (STK17A), Active	S33-10G	Human	GST	Sf9 Cells	Full Length
DRAK2 (STK17B), Active	S34-10G	Human	GST	Sf9 Cells	Full Length
DTYMK, Active	D31-11G	Human	GST	Sf9 Cells	2-end
DYRK1A, Active	D09-10G	Rat	GST	E.coli	Full Length
DYRK1B, Active	D09-10BG	Human	GST	Sf9 Cells	Full-Length
DYRK2, Active	D10-10G	Human	GST	E.coli	Full Length
DYRK3, Active	D11-11G	Human	GST	E.coli	63-end
DYRK4, Active	D12-10G	Human	GST	Sf9 Cells	Full Length
EEF2K, Active	E01-10G	Human	GST	Sf9 Cells	Full Length
EGFR, Active	E10-112G	Human	GST	Sf9 Cells	668-end
EGFR, Active	E10-112H	Human	HIS	Sf9 Cells	668-end
EIF2AK1 (HRI), Active	H07-10G	Human	GST	Sf9 Cells	Full Length
EIF2AK2, Active	P80-11G	Human	GST	Sf9 Cells	252-end
EIF2AK3, Active	E11-11G	Human	GST	E.coli	563-1115
EIF2AK4 (GCN2), Active	E12-11G	Human	GST	Sf9 Cells	192-1024
EPHA1, Active	E13-11G	Mouse	GST	Sf9 Cells	569-end
EPHA1, Active	E13-18G	Human	GST	Sf9 Cells	569-end
EPHA1, Active	E13-18BG	Human	GST	Sf9 Cells	613-892
EPHA2, Active	E14-11G	Human	GST	Sf9 Cells	561-end
EPHA3, Active	E15-11G	Human	GST	Sf9 Cells	571-end
EPHA4, Active	E16-11G	Human	GST	Sf9 Cells	610-887
EPHA4, Active	E16-11BG	Human	GST	Sf9 Cells	570-end
EPHA6, Active	E18-11G	Human	GST	Sf9 Cells	683-end

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
EPHA7, Active	E19-11G	Mouse	GST	Sf9 Cells	580-end
EPHA8 Protein	E20-35G	Human	GST	Sf9 Cells	565-end
EPHB1, Active	E21-11G	Mouse	GST	Sf9 Cells	591-end
EPHB2, Active	E22-11G	Human	GST	Sf9 Cells	570-end
EPHB2, Active	E22-11BG	Human	GST	Sf9 Cells	611-893
EPHB3, Active	E23-11G	Human	GST	Sf9 Cells	585-end
EPHB4, Active	E24-11G	Human	GST	Sf9 Cells	561-end
EPHB4, Active	E24-11H	Human	HIS	Sf9 Cells	561-end
EPHB6 Protein	E26-35G	Human	GST	Sf9 Cells	617-end
ERK1, Active	M29-10U	Human		E.coli	Full Length
ERK1, Active	M29-10G	Human	GST	E.coli	Full Length
ERK1, Unactive	M29-14U	Human	TAG-Free	E.coli	Full Length
ERK1, Unactive	M29-14G	Human	GST	E.coli	Full Length
ERK2, Active	M28-10G	Human	GST	E.coli	Full Length
ERK2, Unactive	M28-14G	Human	GST	E.coli	Full Length
ERK2, Unactive	M28-14U	Human	TAG-Free	E.coli	Full Length
ERK3 (MAPK6), Active	M31-10G	Human	GST	Sf9 Cells	Full Length
ERK4 (MAPK4) Protein	M30-34G	Human	GST	Sf9 Cells	Full Length
ERK5, Active	M32B-10G	Human	GST	Sf9 Cells	Full length
ERK5, Unactive	M32-14G	Human	GST	Sf9 Cells	Full Length
ERK7 (MAPK15), Active	M47-10BG	Mouse	GST	Sf9 Cells	Full Length
ERN1 (IRE1), Active	E31-11G	Human	GST	Sf9 Cells	468-end
ERN1 (IRE1), Active	E31M-11G	Mouse	GST	Sf9 Cells	469-end
ERN2 (IRE2), Active	E32-11G	Human	GST	Sf9 Cells	499-end
FAK, Active	P91-11H	Human	HIS	Sf9 Cells	393-698
FAK, Active	P91-10G	Human	GST	Sf9 Cells	Full Length
FAM20C, Active	F18-11G	Human	GST	Sf9 Cells	23-end
FASTK, Active	F01-10G	Human	GST	Sf9 Cells	Full Length
FER, Active	F16-11G	Mouse	GST	Sf9 Cells	542-end
FES, Active	F03-10G	Human	GST	Sf9 Cells	Full Length
FGFR1 (FLT2), Active	F04-11G	Human	GST	Sf9 Cells	399-822
FGFR2, Active	F05-11G	Human	GST	Sf9 Cells	285-end
FGFR3 (d613-653), Unactive	F06-16HG	Human	GST	Sf9 Cells	d613-653
FGFR3, Active	F06-11G	Human	GST	Sf9 Cells	397-end
FGFR4, Active	F07-11G	Human	GST	Sf9 Cells	460-end
FGR, Active	F10-10G	Human	GST	Sf9 Cells	Full Length
FLT1, Active	F11-11G	Human	GST	Sf9 Cells	784-end
FLT3, Active	F12-11G	Human	GST	Sf9 Cells	571-993
FLT4, Active	F13-11G	Human	GST	Sf9 Cells	800-end
FMS, Active	C74-11G	Human	GST	Sf9 Cells	539-end
FRK, Active	F14-11G	Human	GST	Sf9 Cells	208-end
FYN A, Active	F15-10G	Human	GST	Sf9 Cells	Full Length
FYN C, Unactive	F15-14G	Human	GST	Sf9 Cells	Full Length
GAK, Active	G01-11G	Human	GST	Sf9 Cells	1-396
GCK, Active	M24-10G	Human	GST	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
GLK, Active	M25-11G	Human	GST	Sf9 Cells	1-380
GRK1, Active	R03-10G	Human	GST	Sf9 Cells	Full Length
GRK2, Active	A14-10G	Human	GST	Sf9 Cells	Full Length
GRK3, Active	A15-11G	Human	GST	Sf9 Cells	1-554
GRK4 Protein	G02-31G	Human	GST	Sf9 Cells	140-end
GRK5, Active	G03-10G	Human	GST	Sf9 Cells	Full Length
GRK6, Active	G04-10G	Human	GST	Sf9 Cells	Full Length
GRK7, Active	G05-10G	Human	GST	Sf9 Cells	Full Length
GSK3 alpha, Active	G08-10G	Human	GST	Sf9 Cells	Full Length
GSK3 beta, Active	G09-10G	Human	GST tag	Sf9 Cells	Full Length
GSK3 beta, Active	G09-10H	Human	HIS	Sf9 Cells	Full Length
GSK3 beta, Unactive	G09-14G	Human	GST tag	E.coli	Full Length
GUCY2F Protein	G12-31G	Human	GST	Sf9 Cells	367-end
H11 (HSP22) Protein	H09-34G	Human	GST	Sf9 Cells	Full Length
Haspin (GSG2), Active	G10-11G	Human	GST	Sf9 Cells	470-end
HCK, Active	H02-11G	Human	GST	Sf9 Cells	230-497
HER2, Active	E27-11G	Human	GST	Sf9 Cells	676-end
HER2, Active	E27-11H	Human	HIS	Sf9 Cells	676-end
HER3 Protein	E28-35G	Human	GST	Sf9 Cells	684-1020
HER3 Protein	E28-35H	Human	HIS	Sf9 Cells	667-1020
HER4, Active	E29-11G	Human	GST	Sf9 Cells	682-993
HGK, Active	M26-11G	Human	GST	Sf9 Cells	1-328
HIPK1, Active	H03-11G	Human	GST	Sf9 Cells	156-555
HIPK2, Active	H04-11G	Mouse	GST	Sf9 Cells	153-564
HIPK2, Active	H04-11BG	Human	GST	Sf9 Cells	1-640
HIPK3, Active	H05-13H	Human	HIS	Sf9 Cells	163-562
HIPK4, Active	H06-10G	Human	GST	Sf9 Cells	Full Length
HIPK4, Active	H06-10H	Human	HIS	Sf9 Cells	Full Length
HPK1, Active	M23-11G	Human	GST	Sf9 Cells	1-346
HPK1, Active	M23-11BG	Mouse	GST	Sf9 Cells	1-346
HPK1, Active	M23-10G	Human	GST	Sf9 Cells	Full Length
HUNK, Active	H08-10G	Human	GST	Sf9 Cells	Full Length
ICK, Active	I01-10G	Human	GST	Sf9 Cells	Full Length
IGF1R, Active	I02-11H	Human	HIS	Sf9 Cells	960-end
IKK alpha, Active	C51-10G	Human	GST	Sf9 Cells	Full Length
IKK beta, Active	I03-10BG	Human	GST	Sf9 Cells	1-662
IKK beta, Active	I03-18G	Human	GST	Sf9 Cells	Full Length
IKK beta, Active	I03M-10BG	Mouse	GST	Sf9 Cells	1-662
IKK epsilon, Active	I04-10G	Human	GST	Sf9 Cells	Full Length
IKK gamma Protein	I25-30G	Human	GST	Sf9 Cells	Full Length
InsR, Active	I08-11G	Human	GST	Sf9 Cells	1011-end
IRAK1 (catalytic domain), Active	I09-11G	Human	GST	Sf9 Cells	194-end
IRAK1, Active	I09-10G	Human	GST	Sf9 Cells	Full Length
IRAK2, Active	I10-10BG	Human	GST	Sf9 Cells	Full Length
IRAK3 Protein	I11-34G	Human	GST	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
IRAK4 (catalytic domain), Active	I12-11G	Human	GST	Sf9 Cells	104 a.a. end
IRAK4, Active	I12-10G	Human	GST	Sf9 Cells	Full Length
IRR, Active	I07-11G	Human	GST	Sf9 Cells	945-end
ITK, Active	I13-11G	Human	GST	Sf9 Cells	352-end
ITK, Active	I13-10G	Human	GST	Sf9 Cells	Full Length
JAK1, Active	J01-11G	Human	GST	Sf9 Cells	438-end
JAK2, Active	J02-11G	Human	GST	Sf9 Cells	804-end
JAK2, Active	J02M-11G	Mouse	GST	Sf9 Cells	808-end
JAK3 (JH2 domain) Protein	J03-35G	Human	GST	Sf9 Cells	551-790
JAK3, Active	J03-11G	Human	GST	Sf9 Cells	781-end
JNK1, Active	M33-10G	Mouse	GST	Sf9 Cells	Full Length
JNK1, Unactive	M33-14G	Mouse	GST	E.coli	Full Length
JNK2, Active	M34-10BG	Human	GST	Sf9 Cells	Full Length
JNK2, Unactive	M34-14G	Rat	GST	E.coli	Full Length
JNK3, Active	M35-10BG	Human	GST	Sf9 Cells	Full Length
JNK3, Unactive	M35-14G	Human	GST	E.coli	Full Length
KDR, Active	K01-11G	Human	GST	Sf9 Cells	789-end
KHS1, Active	M27-10G	Human	GST	Sf9 Cells	Full Length
KSR1, Unactive	K07-15BG	Human	GST	Sf9 Cells	335-end
KSR2, Unactive	K08-15G	Human	GST	Sf9 Cells	554-end
LATS1, Active	L01-11G	Human	GST	Sf9 Cells	589-end
LATS2, Active	L02-11G	Human	GST	Sf9 Cells	480-1088
LATS2, Unactive	L02-35G	Human	GST	Sf9 Cells	480-1088
LCK, Active	L03-10G	Human	GST	Sf9 Cells	Full Length
LIMK1, Active	L04-11G	Human	GST	Sf9 Cells	285-638
LIMK1, Unactive	L04-15G	Human	GST	Sf9 Cells	285-638
LIMK1, Unactive	L04-15H	Human	HIS	Sf9 Cells	285-638
LIMK2, Active	L05-10G	Human	GST	Sf9 Cells	Full Length
LKB1(STK11) Protein	L15-30G	Human	GST	Sf9 Cells	Full Length
LMTK2 Protein	L07-35G	Human	GST	Sf9 Cells	66-583
LMTK3 Protein	L08-35G	Mouse	GST	Sf9 Cells	63-484
LOK, Active	S29-11G	Human	GST	Sf9 Cells	1-348
LOK, Active	S29-10G	Human	GST	Sf9 Cells	Full Length
LRRK1, Active	L09-11G	Human	GST	Sf9 Cells	622-1610
LRRK2, Active	L10-11G	Human	GST	Sf9 Cells	968-end
LYN A, Active	L13-18G	Human	GST	Sf9 Cells	Full Length
LYN B, Active	L13-10G	Human	GST	Sf9 Cells	Full Length
LZK, Active	M21-11G	Human	GST	Sf9 Cells	1-427
MAK, Active	M01-10G	Human	GST	Sf9 Cells	Full Length
MAPKAPK2, Active	M40-11G	Human	GST	Sf9 Cells	46-end
MAPKAPK2, Active	M40-11H	Human	HIS	Sf9 Cells	46-end
MAPKAPK2, Unactive	M40-14G	Human	GST	E.coli	46-end
MAPKAPK3, Active	M41-10G	Human	GST	Sf9 Cells	Full Length
MAPKAPK3, Unactive	M41-14G	Human	GST	E.coli	Full Length
MAPKAPK5, Active	M42-10G	Human	GST	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
MAPKAPK5, Unactive	M42-14G	Human	GST	Sf9 Cells	Full Length
MAPKAPK5, Unactive	M42-14BH	Human	HIS	E.coli	Full length
MARK1, Active	M43-10G	Human	GST	Sf9 Cells	Full Length
MARK2, Active	M44-10G	Human	GST	Sf9 Cells	Full Length
MARK3, Active	M45-10G	Human	GST	Sf9 Cells	Full Length
MARK4, Active	M46-10G	Human	GST	Sf9 Cells	Full Length
MAST2, Active	M69-11G	Human	GST	Sf9 Cells	Full Length
MAST3, Active	M70-11G	Human	GST	Sf9 Cells	1-723a.a.
MASTL Protein	M72-30G	Human	GST	Sf9 Cells	Full Length
MASTL, Active	M72-10G	Human	GST	Sf9 Cells	Full Length
MATK (CTK), Active	M49-10G	Human	GST	Sf9 Cells	Full Length
MEK1 (EE), Active	M02-12G	Mouse	GST	E.coli	Full Length
MEK1, Active	M02-10G	Human	GST	Sf9 Cells	Full Length
MEK1, Unactive	M02-14H	Human	HIS	Sf9 Cells	Full Length
MEK1, Unactive	M02-14BG	Human	GST	Sf9 Cells	Full Length
MEK2, Active	M03-10G	Human	GST	Sf9 Cells	Full Length
MEK2, Unactive	M03-14G	Human	GST	Sf9 Cells	Full Length
MEK3 (MAP2K3), Active	M04-10G	Human	GST	Sf9 Cells	Full Length
MEK3 (MAP2K3), Unactive	M04-14G	Human	GST	Sf9 Cells	Full Length
MEK4 (MAP2K4), Active	M05-10G	Human	GST	Sf9 Cells	Full Length
MEK4 (MAP2K4), Unactive	M05-15G	Human	GST	E.coli	33-end
MEK5, Active	M06-10G	Human	GST	Sf9 Cells	Full Length
MEK6 (MAP2K6), Unactive	M07-14G	Human	GST	Sf9 Cells	Full Length
MEK6 (MAP2K6), Unactive	M07-14H	Human	HIS	E.coli	Full Length
MEK6, Active	M07-10G	Human	GST	Sf9 Cells	Full Length
MEK7, Active	M08-10G	Human	GST	Sf9 Cells	Full Length
MEKK1, Active	M09-11G	Human	GST	Sf9 Cells	900-1748
MEKK2, Active	M10-10G	Human	GST	Sf9 Cells	Full Length
MEKK3, Active	M11-10G	Human	GST	Sf9 Cells	Full Length
MEKK4, Active	M12-11G	Human	GST	Sf9 Cells	1203-end
MEKK6 (ASK2), Active	M14-11G	Human	GST tag	Sf9 Cells	620-950
MEKK6 (ASK2), Active	M14-10G	Human		Sf9 Cells	Full Length
MELK, Active	M50-11G	Human	GST	Sf9 Cells	1-340
MELK, Active	M50-18G	Human	GST	Sf9 Cells	1-550
MET, Active	M52-11G	Rat	GST	Sf9 Cells	958-end
MET, Active	M52-18G	Human	GST	Sf9 Cells	956-end
MET, Active	M52-18H	Human	HIS	Sf9 Cells	956-end
MINK1, Active	M53-11G	Human	GST	Sf9 Cells	1-320
MKK4, Active	M05-10H	Human	HIS	Sf9 Cells	Full Length
MLCK, Active	M62-11G	Human	GST	Sf9 Cells	1425-1776
MLK1, Active	M17-11G	Human	GST	Sf9 Cells	1-433
MLK2, Active	M18-11G	Human	GST	Sf9 Cells	1-446
MLK3, Active	M19-11G	Human	GST	Sf9 Cells	1-488
MLK4, Active	M48-11G	Human	GST	Sf9 Cells	114-420
MNK1, Active	M54-10G	Human	GST	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
MNK2, Active	M55-10G	Human	GST	Sf9 Cells	Full Length
MOS Protein	M57-34G	Human	GST	Sf9 Cells	Full Length
MRCK alpha, Active	C27-11G	Human	GST	Sf9 Cells	1-473
MRCK beta, Active	C28-11G	Human	GST	Sf9 Cells	1-473
MSK1, Active	R19-10G	Human	GST	Sf9 Cells	Full Length
MSK2, Active	R18-10G	Mouse	GST	Sf9 Cells	Full Length
MSSK1, Active	S41-10G	Human	GST	Sf9 Cells	Full Length
MST1, Active	S25-10G	Human	GST	Sf9 Cells	Full Length
MST3, Active	S42-11G	Human	GST	Sf9 Cells	1-311
MST4, Active	M59-10G	Human	GST	Sf9 Cells	Full Length
MUSK, Active	M60-11G	Human	GST	Sf9 Cells	519-end
MYLK2, Active	M63-10H	Human	HIS	Sf9 Cells	Full Length
MYLK3, Active	M64-10G	Human	GST	Sf9 Cells	Full Length
MYLK4, Active	M74-10G	Human	GST	Sf9 Cells	Full Length
MYO3 beta, Active	M66-11G	Human	GST	Sf9 Cells	1-326
MYO3alpha, Active	M65-11G	Human	GST	Sf9 Cells	1-434
MYT1, Active	M67-10G	Human	GST	Sf9 Cells	Full Length
MYT1, Active	M67-10H	Human	HIS	Sf9 Cells	Full Length
NDR, Active	S49-10G	Human	GST	Sf9 Cells	Full Length
NDR, Unactive	S49-14G	Human	GST	Sf9 Cells	Full Length
NDR2, Active	S50-10G	Human	GST	Sf9 Cells	Full Length
NEK1, Active	N02-11G	Mouse	GST	Sf9 Cells	1-495
NEK1, Active	N02-18G	Human	GST	Sf9 Cells	1-477
NEK10, Active	N12-11G	Human	GST	Sf9 Cells	266-895
NEK11, Active	N13-10G	Human	GST	Sf9 Cells	Full Length
NEK2, Active	N03-10G	Human	GST	Sf9 Cells	Full Length
NEK3, Active	N05-10G	Human	GST	Sf9 Cells	Full Length
NEK4, Active	N06-10G	Human	GST	Sf9 Cells	Full Length
NEK5, Active	N07-11G	Human	GST	Sf9 Cells	1-508
NEK6, Active	N08-10G	Human	GST	Sf9 Cells	Full Length
NEK7, Active	N09-10G	Human	GST	Sf9 Cells	Full Length
NEK8, Active	N10-11G	Human	GST	Sf9 Cells	1-322
NEK9, Active	N11-11G	Human	GST	Sf9 Cells	1-979; deleted fragment of 346-731
NIK, Active	M22-11G	Human	GST	Sf9 Cells	325-end
NIM1, Active	N14-10G	Human	GST	Sf9 Cells	Full Length
NLK, Active	N15-10G	Human	GST	Sf9 Cells	Full Length
NUAK1, Active	N19-10G	Human	GST	Sf9 Cells	Full Length
NUAK2, Active	N20-10G	Human	GST	Sf9 Cells	Full Length
OSR1 Protein	O09-34G	Human	GST	Sf9 Cells	Full Length
p38 alpha, Active	M39-10BG	Human	GST	Sf9 Cells	Full Length
p38 alpha, Unactive	M39-14G	Human	GST	E.coli	Full Length
p38 alpha, Unactive	M39-14U	Human	TAG-Free	E.coli	Full Length
p38 beta, Active	M36-10BG	Human	GST	Sf9 Cells	Full Length
p38 beta, Unactive	M36-14G	Human	GST	E.coli	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
p38 delta, Active	M38-10BG	Human	GST	Sf9 Cells	Full Length
p38 delta, Unactive	M38-14G	Human	GST	Sf9 Cells	Full Length
p38 gamma, Active	M37-10BG	Human	GST	Sf9 Cells	Full Length
p38 gamma, Unactive	M37-14G	Human	GST	Sf9 Cells	Full Length
p70S6K, Active	R21-10H	Human	HIS	Sf9 Cells	Full Length
p70S6K, Unactive	R21-14H	Human	HIS	Sf9 Cells	Full Length
p70S6Kb, Active	R22-10G	Human	GST	Sf9 Cells	Full Length
p70S6Kb, Unactive	R22-14G	Human	GST	Sf9 Cells	Full Length
p70S6KL1 Protein	R25-34G	Human	GST	Sf9 Cells	Full Length
PAK1, Active	P02B-10BG	Human	GST tag	Sf9 Cells	Full length
PAK1, Unactive	P02-17G	Human	GST	Sf9 Cells	Full Length
PAK1/CDC42, Active	P02-10G	Human		Sf9 Cells	Full Length
PAK2 Protein	P03-34G	Human	GST	Sf9 Cells	Full Length
PAK3, Active	P04-10G	Mouse	GST	Sf9 Cells	Full Length
PAK3, Active	P04-18G	Human	GST	Sf9 Cells	Full Length
PAK4, Active	P05-10BG	Human	GST	Sf9 Cells	Full Length
PAK6, Active	P06-10G	Human	GST	Sf9 Cells	Full Length
PAK7, Active	P07-10G	Human	GST	Sf9 Cells	Full Length
PASK, Active	P08-11G	Human	GST	Sf9 Cells	981-end
PCTK1 (CDK16)/CyclinY, Active	P09-11G	Human	GST	Sf9 Cells	PCTK1 (CDK16) (107-end) and Cyclin Y (2-end)
PCTK1 Protein	P09-35G	Human	GST	Sf9 Cells	107-end
PCTK2 (CDK17)/CyclinY, Active	P10-10G	Human	GST	Sf9 Cells	PCTK2 (CDK17) (Full Length) and Cyclin Y (2-end)
PCTK2 Protein	P10-34G	Human	GST	Sf9 Cells	Full Length
PCTK3 (CDK18)/CyclinY, Active	P11-10G	Human	GST	Sf9 Cells	PCTK3 (CDK18)(Full Length) and Cyclin Y (2-end)
PCTK3 Protein	P11-34G	Human	GST	Sf9 Cells	Full Length
PDGFR alpha, Active	P12-11G	Mouse	GST	Sf9 Cells	550-end
PDGFR alpha, Active	P12-18G	Human	GST	Sf9 Cells	550-end
PDGFR beta, Active	P13-11G	Human	GST	Sf9 Cells	557-end
PDHK1, Active	P57-10G	Human	GST	Sf9 Cells	Full Length
PDHK2, Active	P58-10G	Human	GST	Sf9 Cells	Full Length
PDHK3, Active	P59-10G	Human	GST	Sf9 Cells	Full Length
PDHK4, Active	P60-10G	Human	GST	Sf9 Cells	Full Length
PDK1, Active	P14-10H	Human	HIS	Sf9 Cells	Full Length
PEAK1, Unactive	P18-14G	Human	GST	Sf9 Cells	1127-end
PFTK1 (CDK14) Protein	P15-30G	Human	GST	Sf9 Cells	Full Length
PFTK1 (CDK14)/CyclinY, Active	P15-10G	Human	GST	Sf9 Cells	PFTK1 (CDK14) Full Length and Cyclin Y (2-end)
PHKG1, Active	P19-10G	Human	GST	Sf9 Cells	Full Length
PHKG2, Active	P20-10G	Human	GST	Sf9 Cells	Full Length
PIM1, Active	P35-10G	Human	GST	Sf9 Cells	Full Length
PIM2, Active	P36-10G	Human	GST	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
PIM3, Active	P37-10G	Mouse	GST	Sf9 Cells	Full Length
PIM3, Active	P37-10BG	Human	GST	Sf9 Cells	Full Length
PINK1 Protein	P33-35G	Human	GST	Sf9 Cells	114-end
PIP5K3 Protein	P17-31G	Human	GST	Sf9 Cells	1493-end
PKAc alpha, Active	P51-10G	Human	GST	Sf9 Cells	Full Length
PKAc beta, Active	P52-10G	Human	GST	Sf9 Cells	Full Length
PKAc gamma, Active	P53-10G	Human	GST	Sf9 Cells	Full Length
PKC alpha, Active	P61-10G	Xenopus	GST	Sf9 Cells	Full Length
PKC alpha, Active	P61-18G	Human	GST	Sf9 Cells	Full Length
PKC beta I, Active	P62-10G	Human	GST	Sf9 Cells	Full Length
PKC beta II, Active	P63-10G	Human	GST	Sf9 Cells	Full Length
PKC delta, Active	P64-10G	Human	GST	Sf9 Cells	Full Length
PKC epsilon, Active	P65-10G	Human	GST	Sf9 Cells	Full Length
PKC eta, Active	P67-10G	Human	GST	Sf9 Cells	Full Length
PKC gamma, Active	P66-10G	Human	GST	Sf9 Cells	Full Length
PKC iota, Active	P68-10G	Human	GST	Sf9 Cells	Full Length
PKC mu, Active	P72-10G	Human	GST	Sf9 Cells	Full Length
PKC nu, Active	P73-10G	Human	GST	Sf9 Cells	Full Length
PKC theta, Active	P74-10G	Human	GST	Sf9 Cells	Full Length
PKC zeta, Active	P75-10G	Human	GST	Sf9 Cells	Full Length
PKD2 (PRKD2), Active	P76-10G	Human	GST	Sf9 Cells	Full Length
PKN1/PRK1, Active	P70-11G	Human	GST	Sf9 Cells	289-end
PKN2/PRK2, Active	P71-10G	Human	GST	Sf9 Cells	Full Length
PKN3/PRK3, Active	P69-10G	Human	GST	Sf9 Cells	Full Length
PLK1, Active	P41-10H	Human	HIS	Sf9 Cells	Full Length
PLK2, Active	P42-10G	Human	GST	Sf9 Cells	Full Length
PLK3, Active	P43-13G	Human	GST	Sf9 Cells	57-340
PLK3, Active	P43-13H	Human	HIS	Sf9 Cells	57-340
PLK4, Active	P44-11G	Human	GST	Sf9 Cells	Full Length
PRKG1, Active	P78-10BG	Human	GST	Sf9 Cells	Full Length
PRKG1, Active	P78-11G	Human	GST	Sf9 Cells	324-end
PRKG2, Active	P79-10G	Human	GST	Sf9 Cells	Full Length
PRKX, Active	P81-10G	Human	GST	Sf9 Cells	Full Length
PRKY Protein	P82-35G	Human	GST	Sf9 Cells	40-end
PSKH1 Protein	P89-34G	Human	GST	Sf9 Cells	2-end
PTK7 Protein	P95-35G	Human	GST	Sf9 Cells	726-end
PTK9 Protein	P97-34G	Human	GST	Sf9 Cells	2-end
PTK9L (A6R) Protein	P98-30G	Human	GST	Sf9 Cells	Full Length
PXK Protein	P99-35G	Human	GST	Sf9 Cells	Full Length
PYK2, Active	P92-11H	Human	HIS	Sf9 Cells	360-690
QIK, Active	S15-10G	Human	GST	Sf9 Cells	Full Length
RAF1, Active	R01-10G	Human	GST	Sf9 Cells	306-end
RET, Active	R02-11G	Human	GST	Sf9 Cells	658-end
RET, Active	R02-11H	Human	HIS	Sf9 Cells	658-end
RIOK1 Protein	R04-34G	Human	GST	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
RIOK2 Protein	R05-34G	Human	GST	Sf9 Cells	Full Length
RIOK3 Protein	R06-34G	Human	GST	Sf9 Cells	Full Length
RIPK1, Active	R07-10G	Human	GST	Sf9 Cells	Full Length
RIPK1, Active	R07-11G	Human	GST	Sf9 Cells	1-327
RIPK1, Active	R07M-11G	Mouse	GST	Sf9 Cells	1-327
RIPK1, Active	R07C-11G	Cynomolgus Monkey	GST	Sf9 Cells	1-326
RIPK1, Active	R07D-112G	Dog	GST	Sf9 Cells	1-373
RIPK1, Active	R07C-112G	Monkey	GST	Sf9 Cells	1-373
RIPK1, Active	R07D-11G	Dog	GST	Sf9 Cells	1-327
RIPK2, Active	R08-11G	Human	GST	Sf9 Cells	1-299
RIPK3, Active	R09-10G	Human	GST	Sf9 Cells	Full Length
RIPK4 Protein	R26-35G	Human	GST	Sf9 Cells	15-426
RIPK5, Active	R27-10G	Human	GST	Sf9 Cells	Full Length
ROCK1, Active	R10-11G	Human	GST	Sf9 Cells	17-535
ROCK2, Active	R11-11G	Human	GST	Sf9 Cells	5-554
ROCK2, Active	R11-11H	Human	HIS	Sf9 Cells	5-554
RON, Active	M58-11G	Human	GST	Sf9 Cells	983-end
ROR1, Active	R12-11G	Human	GST	Sf9 Cells	429-end
ROR2, Active	R13-11G	Human	GST	Sf9 Cells	427-end
ROS1, Active	R14-11G	Human	GST	Sf9 Cells	1883-end
ROS1, Active	R14-11BG	Human	GST	Sf9 Cells	1881-end
RSK1, Active	R15-10G	Human	GST	Sf9 Cells	Full Length
RSK2, Active	R17-10G	Human	GST	Sf9 Cells	Full Length
RSK3, Active	R16-10G	Human	GST	Sf9 Cells	Full Length
RSK4, Active	R20-10G	Human	GST	Sf9 Cells	Full Length
RYK Protein	R28-31G	Human	GST	Sf9 Cells	255-end
SBK1, Active	S56-10G	Human	GST	Sf9 Cells	Full Length
SGK1, Active	S06-11G	Human	GST	Sf9 Cells	60-end
SGK1, Active	S06-10G	Human	GST	Sf9 Cells	Full Length
SGK1, Active	S06-11H	Human	HIS	Sf9 Cells	60-end
SGK1, Unactive	S06-15G	Human	GST	Sf9 Cells	60-end
SGK1, Unactive	S06-14G	Human	GST	Sf9 Cells	Full Length
SGK2, Active	S07-10G	Human	GST	Sf9 Cells	Full Length
SGK2, Unactive	S07-14G	Human	GST	Sf9 Cells	Full Length
SGK3, Active	S08-10G	Human	GST	Sf9 Cells	Full Length
SGK3, Unactive	S08-14G	Human	GST	Sf9 Cells	Full Length
SGK494 Protein	S57-34G	Human	GST	Sf9 Cells	Full-Length
SIK, Active	S14-11G	Human	GST	Sf9 Cells	1-303
SIK, Active	S14-11BH	Human	HIS	Sf9 Cells	1-303
SIK3, Active	S12-11G	Human	GST	Sf9 Cells	1-307
SLK, Active	S11-10G	Human	GST	Sf9 Cells	Full Length
SNRK, Active	S16-10G	Human	GST	Sf9 Cells	Full Length
SOK1, Active	S43-10G	Human	GST	Sf9 Cells	Full Length
SOK1, Unactive	S43-14G	Human	GST	Sf9 Cells	Full Length

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
SRC, Active	S19-10G	Rous Sarcoma Virus	GST	Sf9 Cells	Full Length
SRC, Active	S19-18G	Human	GST	Sf9 Cells	Full Length
SRMS, Active	S20-11G	Human	GST	Sf9 Cells	215-end
SRPK1, Active	S21-10G	Human	GST	Sf9 Cells	Full Length
SRPK2, Active	S22-10G	Human	GST	Sf9 Cells	Full Length
STK16 Protein	S31-34G	Human	GST	Sf9 Cells	Full Length
STK19, Active	S35-10G	Human	GST	Sf9 Cells	Full Length
STK21 (CIT), Active	C52-11G	Human	GST	Sf9 Cells	1-449
STK3, Active	S24-10G	Human	GST	Sf9 Cells	Full Length
STK31 Protein	S45-35G	Human	GST	Sf9 Cells	496-end
STK32A (YANK1) Protein	Y03-34G	Mouse	GST	Sf9 Cells	Full Length
STK32B (YANK2), Active	Y04-10G	Human	GST	Sf9 Cells	Full Length
STK32C (YANK3), Active	Y05-11G	Human	GST	Sf9 Cells	66-end
STK33, Active	S46-10G	Human	GST	Sf9 Cells	Full Length
STK35 (CLIK1) Protein	S47-35G	Human	GST	Sf9 Cells	134-end
STK36, Active	S48-11G	Human	GST	Sf9 Cells	1-457
STK39 (STLK3), Active	S51-10H	Human	HIS	Sf9 Cells	Full Length
STK40 (SHIK) Protein	S54-34G	Human	GST	Sf9 Cells	Full Length
STK9 Protein	C46-31G	Human	GST	Sf9 Cells	1-498
STYK1(NOK) Protein	S53-35G	Human	GST	Sf9 Cells	58-end
SYK, Active	S52-10G	Human	GST	Sf9 Cells	Full Length
TAK1-TAB1, Active	M15-13G	Human	GST	Sf9 Cells	TAK1 (1-303) and TAB1 (437-end)), linked by a small peptide (DFGGGGG)
TAOK1, Active	T24-11G	Human	GST	Sf9 Cells	1-314
TAOK2, Active	T25-11G	Human	GST	Sf9 Cells	1-314
TAOK3, Active	T26-11G	Human	GST	Sf9 Cells	1-463
TBCKL Protein	T35-31G	Human	GST	Sf9 Cells	Full Length
TBK1, Active	T02-10G	Human	GST	Sf9 Cells	Full Length
TBK1, Active	T02M-10G	Mouse	GST	Sf9 Cells	Full Length
TEC, Active	T03-10G	Human	GST	Sf9 Cells	Full Length
TESK1 Protein	T05-34G	Human	GST	Sf9 Cells	Full Length
TESK2, Active	T06-10G	Human	GST	Sf9 Cells	Full Length
TGFBR1 (ALK5), Active	T07-11G	Human	GST	Sf9 Cells	80-end
TGFBR1 (ALK5), Active	T07-11BG	Mouse	GST	Sf9 Cells	148-end
TGFBR1, Unactive	T07-35G	Human	GST	Sf9 Cells	80-end
TGFBR2, Active	T08-11G	Human	GST	Sf9 Cells	190-end
TGFBR2, Active	T08-11BG	Mouse	GST	Sf9 Cells	215-end
TIE2, Active	T04-11G	Human	GST	Sf9 Cells	771-end
TLK1, Active	T10-10G	Human	GST	Sf9 Cells	Full Length
TLK2, Active	T11-11G	Human	GST	Sf9 Cells	388-end
TNIK, Active	T27-11G	Human	GST	Sf9 Cells	1-367
TNK1, Active	T12-11G	Human	GST	Sf9 Cells	1-510
TNNI3K, Active	T13-11G	Human	GST	Sf9 Cells	300-end

# Protein Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
TOPK, Active	T14-10G	Human	GST	Sf9 Cells	Full Length
TOPK, Unactive	T14-14G	Human	GST	Sf9 Cells	Full Length
TP53RK Protein	T29-34G	Human	GST	Sf9 Cells	Full Length
TRIB1 Protein	T30-30G	Human	GST	Sf9 Cells	Full Length
TRIB2 Protein	T31-30G	Human	GST	Sf9 Cells	Full Length
TRIB3 Protein	T32-30G	Human	GST	Sf9 Cells	Full Length
TRIO, Active	T16-11G	Human	GST	Sf9 Cells	2627-end
TRKA, Active	N16-11G	Human	GST	Sf9 Cells	440-end
TRKB, Active	N17-11G	Human	GST	Sf9 Cells	455-end
TRKC, Active	N18-11H	Human	HIS	Sf9 Cells	507-end
TSSK1B, Active	S40-10BG	Human	GST	Sf9 Cells	Full Length
TSSK2, Active	S38-10G	Human	GST	Sf9 Cells	Full Length
TSSK3 Protein	S39-34G	Human	GST	Sf9 Cells	2-end
TSSK6 (SSTK) Protein	S23-34G	Human	GST	Sf9 Cells	Full Length
TTBK1, Active	T17-11G	Human	GST	Sf9 Cells	1-479
TTBK2, Active	T18-11G	Mouse	GST	Sf9 Cells	70-538
TTK, Active	T20-10G	Human	GST	Sf9 Cells	Full Length
TTK, Active	T20M-10G	Mouse	GST	Sf9 Cells	Full Length
TXK, Active	T19-11G	Human	GST	Sf9 Cells	239-end
TYK2 (JTK1), Active	T21-11G	Human	GST	Sf9 Cells	442-end
TYR03, Active	T22-11G	Human	GST	Sf9 Cells	455-end
UHMK1 (KIS) Protein	K05-30G	Human	GST	Sf9 Cells	Full Length
ULK1, Active	U01-11G	Human	GST	Sf9 Cells	1-649
ULK2, Active	U02-11G	Human	GST	Sf9 Cells	1-478
ULK3, Active	U03-10H	Human	HIS	Sf9 Cells	Full Length
ULK4 Protein	U04-35G	Human	GST	Sf9 Cells	1-616
VLK Protein	V05-31G	Human	GST	Sf9 Cells	129-end
VRK1, Active	V01-10G	Human	GST	Sf9 Cells	Full Length
VRK2, Active	V02-11G	Human	GST	Sf9 Cells	1-375
VRK3 Protein	V03-30G	Human	GST	Sf9 Cells	Full Length
WEE1, Active	W01-11G	Human	GST	Sf9 Cells	Full Length
WEE2, Active	W06-10G	Mouse	GST	Sf9 Cells	Full Length
WNK1, Active	W02-11G	Human	GST	Sf9 Cells	181-507
WNK2, Active	W03-11G	Human	GST	Sf9 Cells	1-489
WNK3, Active	W04-11G	Human	GST	Sf9 Cells	1-571
WNK4, Active	W05-11G	Mouse	GST	Sf9 Cells	1-441
YES1, Active	Y01-10G	Human	GST	Sf9 Cells	Full Length
YSK4 (MAP3K19), Active	Y06-11G	Human	GST	Sf9 Cells	941-end
ZAK, Active	Z01-10G	Human	GST	Sf9 Cells	Full Length
ZAP70, Active	Z02-10G	Human	GST	Sf9 Cells	Full Length

# Lipid Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
DGKA, Active	D21-10BG	Human	GST	Sf9 Cells	Full Length
DGKA, Active	D21-10BH	Human	HIS	Sf9 Cells	Full Length
DGKA, Active	D21-10CH	Rat	HIS	Sf9 Cells	Full Length
DGKA, Active	D21-10DH	Dog	HIS	Sf9 Cells	Full Length
DGKA, Active	D21-10EH	Monkey	HIS	Sf9 Cells	Full Length
DGKA, Active	D21-10CG	Mouse	GST	Sf9 Cells	Full Length
DGKA, Active	D21-10FH	Mouse	HIS	Sf9 Cells	Full Length
DGKB (11-E), Active	D22-11H	Human	HIS	Sf9 Cells	11-end
DGKB, Active	D22-10G	Human	GST	Sf9 Cells	Full Length
DGKB, Active	D22-10H	Human	HIS	Sf9 Cells	Full Length
DGKD, Active	D23-10G	Human	GST	Sf9 Cells	Full Length
DGKD, Active	D23-10H	Human	HIS	Sf9 Cells	Full Length
DGKE, Active	D24-10G	Human	GST	Sf9 Cells	Full Length
DGKE, Active	D24-10H	Human	HIS	Sf9 Cells	Full Length
DGKG, Active	D25-10G	Human	GST	Sf9 Cells	Full Length
DGKG, Active	D25-10H	Human	HIS	Sf9 Cells	Full Length
DGKG, Active	D25-10BH	Human	HIS	Sf9 Cells	Full Length
DGKG, Active	D25-10BG	Human	GST	Sf9 Cells	Full Length
DGKG, Active	D25-11H	Human	HIS	Sf9 Cells	10-end
DGKH, Active	D26-10G	Human	GST	Sf9 Cells	Full Length
DGKH, Active	D26-10H	Human	HIS	Sf9 Cells	Full Length
DGKI, Active	D27-10G	Human	GST	Sf9 Cells	Full Length
DGKI, Active	D27-10H	Human	HIS	Sf9 Cells	Full Length
DGKK, Active	D28-10G	Human	GST	Sf9 Cells	Full Length
DGKK, Active	D28-10H	Human	HIS	Sf9 Cells	Full Length
DGKQ, Active	D29-10G	Human	GST	Sf9 Cells	Full Length
DGKZ, Active	D30-10G	Human	GST	Sf9 Cells	Full Length
DGKZ, Active	D30-10H	Human	HIS	Sf9 Cells	Full Length
DGKZ, Active	D30M-10H	Mouse	HIS	Sf9 Cells	Full Length
DGKZ, Active	D30A-10H	Rat	HIS	Sf9 Cells	Full Length
PI3K (p110 alpha/p55 gamma), Active	P27-10CG	Human	GST	Sf9 Cells	Full Length
PI3K (p110 alpha/p55 gamma), Active	P27-102CH	Human	HIS	Sf9 Cells	Full Length
PI3K (p110 alpha/p85 alpha), Active	P27-102G	Human	GST	Sf9 Cells	Full Length
PI3K (p110 alpha/p85 alpha), Active	P27-102H	Human	HIS	Sf9 Cells	Full Length
PI3K (p110 alpha/p85 alpha), Active	P27-182H	Mouse	HIS	Sf9 Cells	Full Length
PI3K (p110 beta/p85 alpha), Active	P28-102H	Human	HIS	Sf9 Cells	Full Length
PI3K (p110 beta/p85 beta), Active	P28-102BH	Human	HIS	Sf9 Cells	Full Length
PI3K (p110 delta/p85 alpha), Active	P30-102H	Human	HIS	Sf9 Cells	Full Length
PI3K (p120 gamma), Active	P29-102H	Human	HIS	Sf9 Cells	Full Length
PI4K2A, Active	P21-102G	Human	GST	Sf9 Cells	Full Length
PI4K2B, Active	P22-102G	Human	GST	Sf9 Cells	Full Length
PI4KA, Active	P31-102G	Human	GST	Sf9 Cells	Full Length
PI4KB, Active	P32-102G	Human	GST	Sf9 Cells	Full Length
PIK3C2A, Active	P23-112G	Human	GST	Sf9 Cells	299-end
PIK3C2B, Active	P24-112BG	Human	GST	Sf9 Cells	850-end

# Lipid Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
PIK3C2G, Active	P25-112G	Human	GST	Sf9 Cells	468-1203
PIK3C3 (Vps34), Active	P26-102G	Human	GST	Sf9 Cells	Full Length
PIP4K2A, Active	P76-102AG	Human	GST	Sf9 Cells	Full Length
PIP4K2B, Active	P76-102BG	Human	GST	Sf9 Cells	Full Length
PIP4K2C, Active	P76-102CG	Human	GST	Sf9 Cells	Full Length
PIP5K1A, Active	P16-102AG	Human	GST	Sf9 Cells	Full Length
PIP5K1B, Active	P16-102BG	Human	GST	Sf9 Cells	Full Length
PIP5K1C, Active	P16-102CG	Human	GST	Sf9 Cells	Full Length
PIP5K3 (PIKFYVE), Active	P17-112BG	Human	GST	Sf9 Cells	1400-end
SPHK1, Active	S17-102G	Human	GST	Sf9 Cells	Full Length
SPHK2, Active	S18-102G	Human	GST	Sf9 Cells	Full Length

# Carbohydrate Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
PFKFB1, Active	P321-10G	Human	GST	Sf9 Cells	Full length
PFKFB2 Protein	P322-30G	Human	GST	Sf9 Cells	Full Length
PFKFB3, Active	P323-10G	Human	GST	Sf9 Cells	Full length
PFKFB4, Active	P324-11G	Human	GST	Sf9 Cells	Full length
POMK (SGK196) Protein	S55-34G	Human	GST	Sf9 Cells	Full-Length
Uridylate Kinase	UK05-E311H	E.Coli	HIS	E.coli	Full Length

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
ABL1 (E255K), Active	A03-12BG	Human	GST	Sf9 Cells	27-end (E255K)
ABL1 (E255V), Active	A03-12HH	Human	HIS	Sf9 Cells	27-end (E255V)
ABL1 (F317I), Active	A03-12JG	Human	GST	Sf9 Cells	27-end (F317I)
ABL1 (F317L), Active	A03-12IG	Human	GST	Sf9 Cells	27-end (F317L)
ABL1 (H396P), Active	A03-12LG	Human	GST	Sf9 Cells	27-end (H396P)
ABL1 (K290R) Protein	A03-16NG	Human	HIS	Sf9 Cells	27-end (K290R)
ABL1 (M351T), Active	A03-12KG	Human	GST	Sf9 Cells	27-end (M351T)
ABL1 (Q252H), Active	A03-12FG	Human	GST	Sf9 Cells	27-end (Q252H)
ABL1 (T315I), Active	A03-12DG	Human	GST	Sf9 Cells	27-end (T315I)
ABL1 (V299L), Active	A03-12MG	Human	GST	Sf9 Cells	27-end (V299L)
ABL1 (Y245F Y412F), Active	A03-12QH	Human	HIS	Sf9 Cells	27-end (Y245F Y412F)
ABL1 (Y245F), Active	A03-12PH	Human	HIS	Sf9 Cells	27-end (Y245F)
ABL1 (Y253F), Active	A03-12EG	Human	GST	Sf9 Cells	27-end (Y253F)
ABL1 (Y253H), Active	A03-12GH	Human	HIS	Sf9 Cells	27-end (Y253H)
AKT1 (E17K), Active	A16-12G	Human	GST	Sf9 Cells	Full Length (E17K)
AKT1 (E17K), Active	A16-12H	Human	HIS	Sf9 Cells	Full Length (E17K)
AKT2 (E17K), Active	A17-12G	Human	GST	Sf9 Cells	Full Length (E17K)

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
<b>AKT2 (R274H), Active</b>	A17-12BG	Human	GST	Sf9 Cells	Full Length (R274H)
<b>AKT3 (E17K), Active</b>	A18-12G	Human	GST	Sf9 Cells	Full Length (E17K)
<b>AKT3 (G171R), Active</b>	A18-12BG	Human	GST	Sf9 Cells	Full Length (G171R)
<b>ALK (C1156Y), Active</b>	A19-12DG	Human	GST	Sf9 Cells	1060-end (C1156Y)
<b>ALK (F1174L), Active</b>	A19-12EG	Human	GST	Sf9 Cells	1060-end (F1174L)
<b>ALK (F1174S), Active</b>	A19-12FG	Human	GST	Sf9 Cells	1060-end (F1174S)
<b>ALK (G1202R), Active</b>	A19-12HG	Human	GST	Sf9 Cells	1060-end (G1202R)
<b>ALK (G1269A), Active</b>	A19-12JG	Human	GST	Sf9 Cells	1060-end (G1269A)
<b>ALK (G1269S), Active</b>	A19-12KG	Human	GST	Sf9 Cells	1060-end (G1269S)
<b>ALK (I1171N D1203N), Active</b>	A19-12PG	Human	GST	Sf9 Cells	1060-end (I1171N D1203N)
<b>ALK (L1152R), Active</b>	A19-12CG	Human	GST	Sf9 Cells	1060-end (L1152R)
<b>ALK (L1196M G1202R), Active</b>	A19-12NG	Human	GST	Sf9 Cells	1060-end (L1196M G1202R)
<b>ALK (L1196M), Active</b>	A19-12GG	Human	GST	Sf9 Cells	1060-end (L1196M)
<b>ALK (R1275Q), Active</b>	A19-12LG	Human	GST	Sf9 Cells	1060-end (R1275Q)
<b>ALK (S1206R), Active</b>	A19-12IG	Human	GST	Sf9 Cells	1060-end (S1206R)
<b>ALK (T1151M), Active</b>	A19-12BG	Human	GST	Sf9 Cells	1060-end (T1151M)
<b>ALK (T1151_L1152insT), Active</b>	A19-12G	Human	GST	Sf9 Cells	1060-end (T1151_L1152insT)
<b>ALK2 (G328E), Active</b>	A06-12GG	Human	GST	Sf9 Cells	147-end (G328E)
<b>ALK2 (G328V), Active</b>	A06-12HG	Human	GST	Sf9 Cells	147-end (G328V)
<b>ALK2 (G328W), Active</b>	A06-12IG	Human	GST	Sf9 Cells	147-end (G328W)
<b>ALK2 (G356D), Active</b>	A06-12JG	Human	GST	Sf9 Cells	147-end (G356D)
<b>ALK2 (PF197_198L), Active</b>	A06-12FG	Human	GST	Sf9 Cells	147-end (PF197_198L)
<b>ALK2 (Q207D), Active</b>	A06-13CG	Human	GST	Sf9 Cells	147-end (Q207D)
<b>ALK2 (Q207E), Active</b>	A06-13DG	Human	GST	Sf9 Cells	147-end (Q207E)
<b>ALK2 (R202I), Active</b>	A06-12EG	Human	GST	Sf9 Cells	147-end ((R202I)
<b>ALK2 (R206H), Active</b>	A06-12BG	Human	GST	Sf9 Cells	147-end (R206H)
<b>ALK2 (R206H), Active</b>	A06M-12BG	Mouse	GST	Sf9 Cells	147-end (R206H)
<b>ALK2 (R258G), Active</b>	A06-12KG	Human	GST	Sf9 Cells	147-end (R258G)
<b>ALK3 (Q233D), Active</b>	B04-12BG	Human	GST	Sf9 Cells	187-end (Q233D)
<b>ALK3 (R486Q), Active</b>	B04-12CG	Human	GST	Sf9 Cells	187-end (R486Q)
<b>ANXA4-PKN1(Aex2Pex13), Active</b>	P70-19G	Human	GST	Sf9 Cells	1-3a.a., exons1-2 (ANXA4) - 578a.a.-end, exons13-22 (PKN1)
<b>AURORA B (G160L), Active</b>	A31-10BG	Human	GST	Sf9 Cells	Full Length (G160L)
<b>AXL (R499C), Active</b>	A34-12BG	Human	GST	Sf9 Cells	473-end (R499C)
<b>BCR-RET, Active</b>	R02-19EG	Human	GST	Sf9 Cells	1-426 (BCR) - 713-end (RET)
<b>BRAF (d485-489 P490Y), Active</b>	B08-12IG	Human	GST	Sf9 Cells	381-end or exon 9-18 (d485-489 P490Y)
<b>BRAF (D594G) Protein</b>	B08-12MG	Human	GST	Sf9 Cells	381-end (D594G)
<b>BRAF (G464V), Active</b>	B08-12JG	Human	GST	Sf9 Cells	381-end or exon 9-18 (G464V)
<b>BRAF (G469A), Active</b>	B08-12FG	Human	GST	Sf9 Cells	381-end or exon 9-18 (G469A)
<b>BRAF (K601E), Active</b>	B08-12BG	Human	GST	Sf9 Cells	416-end (K601E)
<b>BRAF (L597V), Active</b>	B08-12HG	Human	GST	Sf9 Cells	381-end or exon 9-18 (L597V)

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
<b>BRAF (R506_K507insVLR), Active</b>	B08-13CG	Human	GST	Sf9 Cells	381-end or exon 9-18 (R506_K507insVLR)
<b>BRAF (T599_V600insT), Active</b>	B08-13BG	Human	GST	Sf9 Cells	416-end (T599_V600insT)
<b>BRAF (V600A), Active</b>	B08-12EG	Human	GST	Sf9 Cells	381-end or exon 9-18 (V600A)
<b>BRAF (V600D), Active</b>	B08-12CG	Human	GST	Sf9 Cells	381-end or exon 9-18 (V600D)
<b>BRAF (V600E), Active</b>	B08-12G	Human	GST	Sf9 Cells	416-end (V600E)
<b>BRAF (V600K), Active</b>	B08-12DG	Human	GST	Sf9 Cells	381-end or exon 9-18 (V600K)
<b>BTK (C481F), Active</b>	B10-12FH	Human	HIS	Sf9 Cells	Full Length (C481F)
<b>BTK (C481R), Active</b>	B10-12GH	Human	HIS	Sf9 Cells	Full Length (C481R)
<b>BTK (C481S), Active</b>	B10-12CH	Human	HIS	Sf9 Cells	Full Length (C481S)
<b>BTK (P190K), Active</b>	B10-12BH	Human	HIS	Sf9 Cells	Full Length (P190K)
<b>BTK (T474I), Active</b>	B10-12HH	Human	HIS	Sf9 Cells	Full Length (T474I)
<b>BTK (Y485F), Active</b>	B10-12IH	Human	HIS	Sf9 Cells	Full Length (Y485F)
<b>c-KIT (A829P), Active</b>	C06-12HG	Human	GST	Sf9 Cells	544-end (A829P)
<b>c-KIT (d557-558 V654A), Active</b>	K06-12B2G	Human	GST	Sf9 Cells	544-end (d557-558 V654A)
<b>c-KIT (d557-558 Y823D), Active</b>	K06-12A2G	Human	GST	Sf9 Cells	544-end (d557-558 Y823D)
<b>c-KIT (d557-558), Active</b>	K06-12WG	Human	GST	Sf9 Cells	544-end (d557-558)
<b>c-KIT (d560-576 Y823D), Active</b>	K06-12YG	Human	GST	Sf9 Cells	544-end (d560-576 Y823D)
<b>c-KIT (D816E), Active</b>	C06-12IG	Human	GST	Sf9 Cells	544-end (D816E)
<b>c-KIT (D816F), Active</b>	C06-12JG	Human	GST	Sf9 Cells	544-end (D816F)
<b>c-KIT (D816H), Active</b>	C06-12KG	Human	GST	Sf9 Cells	544-end (D816H)
<b>c-KIT (D816I), Active</b>	K06-12XG	Human	GST	Sf9 Cells	544-end (D816I)
<b>c-KIT (D816V), Active</b>	C06-12LG	Human	GST	Sf9 Cells	544-end (D816V)
<b>c-KIT (D816Y), Active</b>	C06-12MG	Human	GST	Sf9 Cells	544-end (D816Y)
<b>c-KIT (D820E), Active</b>	K06-12BG	Human	GST	Sf9 Cells	544-end (D820E)
<b>c-KIT (D820Y), Active</b>	K06-12CG	Human	GST	Sf9 Cells	544-end (D820Y)
<b>c-KIT (K642E), Active</b>	C06-12NG	Human	GST	Sf9 Cells	544-end (K642E)
<b>c-KIT (L576P), Active</b>	K06-12UG	Human	GST	Sf9 Cells	544-end (L576P)
<b>c-KIT (N822K), Active</b>	K06-12DG	Human	GST	Sf9 Cells	544-end (N822K)
<b>c-KIT (T670E), Active</b>	K06-12EG	Human	GST	Sf9 Cells	544-end (T670E)
<b>c-KIT (T670I), Active</b>	C06-12FG	Human	GST	Sf9 Cells	544-end (T670I)
<b>c-KIT (V559A), Active</b>	C06-12OG	Human	GST	Sf9 Cells	544-end (V559A)
<b>c-KIT (V559D T670I), Active</b>	K06-12QG	Human	GST	Sf9 Cells	544-end (V559D T670I)
<b>c-KIT (V559D V654A), Active</b>	K06-12RG	Human	GST	Sf9 Cells	544-end (V559D V654A)
<b>c-KIT (V559D), Active</b>	C06-12PG	Human	GST	Sf9 Cells	544-end (V559D)
<b>c-KIT (V560G D816V), Active</b>	C06-12TG	Human	GST	Sf9 Cells	544-end (V560G D816V)
<b>c-KIT (V560G N822K), Active</b>	K06-12VG	Human	GST	Sf9 Cells	544-end (V560G N822K)
<b>c-KIT (V560G), Active</b>	C06-12SG	Human	GST	Sf9 Cells	544-end (V560G)
<b>c-KIT (V654A), Active</b>	K06-13G	Human	GST	Sf9 Cells	544-end (V654A)
<b>c-KIT (Y823D), Active</b>	K06-12GG	Human	GST	Sf9 Cells	544-end (Y823D)
<b>c-MER (A708S), Active</b>	M51-12BG	Human	GST	Sf9 Cells	578-872 (A708S)
<b>CK1 epsilon (R178C), Active</b>	C66-12BG	Human	GST	Sf9 Cells	Full Length (R178C)

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
CK1A1 (E42C), Active	C64-12CG	Human	GST	Sf9 Cells	Full length (E42C)
CK1A1 (I35C), Active	C64-12BG	Human	GST	Sf9 Cells	Full length (I35C)
DDR1 (S496A), Active	D05-12G	Human	GST	Sf9 Cells	444-end (S496A)
DDR2 (T654M), Active	D06-12G	Human	GST	Sf9 Cells	467-end (T654M)
DNAJB1-PRKACA, Active	P51-19BG	Human	GST	Sf9 Cells	2-70 (DNAJB1) - 16-351 (PRKACA)
EGFR (A763_Y764insFHEA), Active	E10-132BG	Human	GST	Sf9 Cells	668-end (A763_Y764insFHEA)
EGFR (A763_Y764insFQEA), Active	E10-132UG	Human	GST	Sf9 Cells	668-end (A763_Y764insFQEA)
EGFR (A767-S768insTLA), Active	E10-132CG	Human	GST	Sf9 Cells	668-end (A767-S768insTLA)
EGFR (C775S T790M L858R), Active	E10-122D2G	Human	GST	Sf9 Cells	668-end (C775S T790M L858R)
EGFR (C797A), Active	E10-12RG	Human	GST	Sf9 Cells	695-end (C797A)
EGFR (C797G L858R), Active	E10-122LG	Human	GST	Sf9 Cells	668-end (C797G L858R)
EGFR (C797S L858R), Active	E10-122ZG	Human	GST	Sf9 Cells	668-end (C797S L858R)
EGFR (C797S), Active	E10-122SG	Human	GST	Sf9 Cells	668-end (C797S)
EGFR (d746), Active	E10-12IG	Human	GST	Sf9 Cells	695-end (d746)
EGFR (d746-750 C775S T790M L858R), Active	E10-122E2G	Human	GST	Sf9 Cells	668-end (d746-750 C775S T790M L858R)
EGFR (d746-750 C797A), Active	E10-122C2G	Human	GST	Sf9 Cells	668-end (d746-750 C797A)
EGFR (d746-750 C797G), Active	E10-122NG	Human	GST	Sf9 Cells	668-end (d746-750 C797G)
EGFR (d746-750 C797S), Active	E10-122TG	Human	GST	Sf9 Cells	668-end (d746-750 C797S)
EGFR (D746-750 G724S), Active	E10-122D3G	Human	GST	Sf9 Cells	668-end (D746-750 G724S)
EGFR (d746-750 L718Q T790M), Active	E10-122E3G	Human	GST	Sf9 Cells	668-end (d746-750 L718Q T790M)
EGFR (d746-750 S768I), Active	E10-122B3G	Human	GST	Sf9 Cells	668-end (d746-750 S768I)
EGFR (d746-750 T790M C797G) Protein	E10-13TG	Human	GST	Sf9 Cells	668-end (d746-750 T790M C797G)
EGFR (d746-750 T790M C797S L858R), Active	E10-122YG	Human	GST	Sf9 Cells	668-end (d746-750 T790M C797S L858R)
EGFR (d746-750 T790M C797S), Active	E10-122UG	Human	GST	Sf9 Cells	668-end (d746-750 T790M C797S)
EGFR (d746-750 T790M L792F), Active	E10-122G3G	Human	GST	Sf9 Cells	668-end (d746-750 T790M L792F)
EGFR (d746-750 T790M L792H), Active	E10-122F3G	Human	GST	Sf9 Cells	668-end (d746-750 T790M L792H)
EGFR (d746-750 T790M L798I), Active	E10-122WG	Human	GST	Sf9 Cells	668-end (d746-750 T790M L798I)
EGFR (d746-750 T790M L858R), Active	E10-122XG	Human	GST	Sf9 Cells	668-end (d746-750 T790M L858R)
EGFR (d746-750 T790M), Active	E10-122KG	Human	GST	Sf9 Cells	668-end (d746-750 T790M)
EGFR (d746-750), Active	E10-122JG	Human	GST	Sf9 Cells	668-end (d746-750)
EGFR (d747-749 A750P), Active	E10-12MG	Human	GST	Sf9 Cells	695-end (d747-749 A750P)
EGFR (d747-749), Active	E10-12LG	Human	GST	Sf9 Cells	695-end (d747-749)
EGFR (d747-752 P753S), Active	E10-12NG	Human	GST	Sf9 Cells	695-end (d747-752 P753S)
EGFR (d752-759), Active	E10-12OG	Human	GST	Sf9 Cells	695-end (d752-759)
EGFR (D761Y), Active	E10-12QG	Human	GST	Sf9 Cells	695-end (D761Y)
EGFR (D761Y), Active	E10-122QG	Human	GST	Sf9 Cells	668-end (D761Y)
EGFR (D770GY), Active	E10-13KG	Human	GST	Sf9 Cells	695-end (D770GY)
EGFR (D770GY), Active	E10-132KG	Human	GST	Sf9 Cells	668-end (D770GY)
EGFR (D770_N771insG), Active	E10-132XG	Human	GST	Sf9 Cells	668-end (D770_N771insG)
EGFR (D770_N771insGF) Protein	E10-13IG	Human	GST	Sf9 Cells	695-end (D770_N771insGF)

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
EGFR (D770_N771insNPG T790M), Active	E10-132TG	Human	GST	Sf9 Cells	668-end (D770_N771insNPG T790M)
EGFR (D770_N771insNPG), Active	E10-132GG	Human	GST	Sf9 Cells	668-end (D770_N771insNPG)
EGFR (D770_N771insY), Active	E10-122D4G	Human	GST	Sf9 Cells	669-end (D770_N771insY)
EGFR (G719C), Active	E10-12FG	Human	GST	Sf9 Cells	695-end (G719C)
EGFR (G719D), Active	E10-12GG	Human	GST	Sf9 Cells	695-end (G719D)
EGFR (G719S), Active	E10-12HG	Human	GST	Sf9 Cells	695-end (G719S)
EGFR (H773_V774insAH) Protein	E10-13PG	Human	GST	Sf9 Cells	695-end (H773_V774insAH)
EGFR (H773_V774insH) Protein	E10-13RG	Human	GST	Sf9 Cells	695-end (H773_V774insH)
EGFR (K716A C797S L858R), Active	E10-122S2G	Human	GST	Sf9 Cells	668-end (K716A C797S L858R)
EGFR (K716A T790M C797S L858R), Active	E10-122P2G	Human	GST	Sf9 Cells	668-end (K716A T790M C797S L858R)
EGFR (K716A), Active	E10-122M2G	Human	GST	Sf9 Cells	668-end (K716A)
EGFR (K716Q L718Q), Active	E10-122T2G	Human	GST	Sf9 Cells	668-end (K716Q L718Q)
EGFR (K728A T790M C797S L858R), Active	E10-122Q2G	Human	GST	Sf9 Cells	668-end (K728A T790M C797S L858R)
EGFR (K728A), Active	E10-122N2G	Human	GST	Sf9 Cells	668-end (K728A)
EGFR (L718Q C797S L858R), Active	E10-122A3G	Human	GST	Sf9 Cells	668-end (L718Q C797S L858R)
EGFR (L718Q L858R), Active	E10-122C3G	Human	GST	Sf9 Cells	668-end (L718Q L858R)
EGFR (L718Q T790M L858R), Active	E10-122H2G	Human	GST	Sf9 Cells	668-end (L718Q T790M L858R)
EGFR (L718Q), Active	E10-122G2G	Human	GST	Sf9 Cells	668-end (L718Q)
EGFR (L747S), Active	E10-12PG	Human	GST	Sf9 Cells	695-end (L747S)
EGFR (L747S), Active	E10-122Y2G	Human	GST	Sf9 Cells	668-end (L747S)
EGFR (L792F L858R), Active	E10-122U2G	Human	GST	Sf9 Cells	668-end (L792F L858R)
EGFR (L792F), Active	E10-122I2G	Human	GST	Sf9 Cells	668-end (L792F)
EGFR (L792H C797S L858R), Active	E10-122V2G	Human	GST	Sf9 Cells	668-end (L792H C797S L858R)
EGFR (L792H L858R), Active	E10-122MG	Human	GST	Sf9 Cells	668-end (L792H L858R)
EGFR (L792H), Active	E10-122J2G	Human	GST	Sf9 Cells	668-end (L792H)
EGFR (L858R), Active	E10-122BG	Human	GST	Sf9 Cells	668-end (L858R)
EGFR (L861Q), Active	E10-12CG	Human	GST	Sf9 Cells	695-end (L861Q)
EGFR (N771-P772insH), Active	E10-132WG	Human	GST	Sf9 Cells	668-end (N771-P772insH)
EGFR (N771_P772insN), Active	E10-132YG	Human	GST	Sf9 Cells	668-end (N771_P772insN)
EGFR (P772_H773insNPH), Active	E10-13MG	Human	GST	Sf9 Cells	695-end (P772_H773insNPH)
EGFR (P772_H773insPR), Active	E10-132ZG	Human	GST	Sf9 Cells	668-end (P772_H773insPR)
EGFR (R999A), Active	E10-122O2G	Human	GST	Sf9 Cells	668-end (R999A)
EGFR (S768I), Active	E10-122I3G	Human	GST	Sf9 Cells	668-end (S768I)
EGFR (T790M C797G L858R), Active	E10-122OG	Human	GST	Sf9 Cells	668-end (T790M C797G L858R)
EGFR (T790M C797S L858R R999A), Active	E10-122R2G	Human	GST	Sf9 Cells	668-end (T790M C797S L858R R999A)
EGFR (T790M C797S L858R), Active	E10-122VG	Human	GST	Sf9 Cells	668-end (T790M C797S L858R)
EGFR (T790M C797S), Active	E10-122B2G	Human	GST	Sf9 Cells	668-end (T790M C797S)
EGFR (T790M L792F C797S L858R), Active	E10-122W2G	Human	GST	Sf9 Cells	668-end (T790M L792F C797S L858R)
EGFR (T790M L792F L858R), Active	E10-122K2G	Human	GST	Sf9 Cells	668-end (T790M L792F L858R)

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
EGFR (T790M L792H C797S L858R), Active	E10-122X2G	Human	GST	Sf9 Cells	668-end (T790M L792H C797S L858R)
EGFR (T790M L792H L858R), Active	E10-122L2G	Human	GST	Sf9 Cells	668-end (T790M L792H L858R)
EGFR (T790M L858R), Active	E10-122DG	Human	GST	Sf9 Cells	668-end (T790M L858R)
EGFR (T790M), Active	E10-122EG	Human	GST	Sf9 Cells	668-end (T790M)
EGFR (V769-D770insGE), Active	E10-132VG	Human	GST	Sf9 Cells	668-end (V769-D770insGE)
EGFR (V774_C775insHV) Protein	E10-13SG	Human	GST	Sf9 Cells	668-end (V774_C775insHV)
EGFR (V948R), Active	E10-122H3H	Human	HIS	Sf9 Cells	669-1022 (V948R)
EML4-ALK (Eex13Aex20), Active	A19-19G	Human	GST	Sf9 Cells	EML4 (1-496) - ALK (1058-end)
EML4-ALK (Eex14Aex20), Active	A19-19GG	Human	GST	Sf9 Cells	EML4 (1-534) - ALK (1058-end)
EML4-ALK (Eex20Aex20), Active	A19-19HG	Human	GST	Sf9 Cells	EML4 (1-580) - ALK (1058-end)
EML4-ALK (Eex6Aex20), Active	A19-19FG	Human	GST	Sf9 Cells	EML4 (1-222) - ALK (1058-end)
ERN1 (R727A R728A), Active	E31-12DG	Human		Sf9 Cells	468-end (R727A R728A)
ERN1 (R727A), Active	E31-12BG	Human	GST	Sf9 Cells	468-end (R727A)
ERN1 (R728A), Active	E31-12CG	Human	GST	Sf9 Cells	468-end (R728A)
FAM131B-BRAF (Fex2Bex9), Active	B08-19EG	Human	GST	Sf9 Cells	1-30 exon 2 (FAB131B) - 381-end or exon 9-18 (BRAF)
FGFR1 (V561M), Active	F04-13G	Human	GST	Sf9 Cells	399-822 (V561M)
FGFR1 (W666R), Active	F04-12G	Human	GST	Sf9 Cells	399-end (W666R)
FGFR10P-FGFR1, Active	F04-19BG	Human	GST	Sf9 Cells	1-173 (FGFR10P) 429-end (FGFR1)
FGFR2 (C491A V564I), Active	F05-12PG	Human	GST	Sf9 Cells	285-end (C491A V564I)
FGFR2 (C491A V564L), Active	F05-12QG	Human	GST	Sf9 Cells	285-end (C491A V564L)
FGFR2 (C491A), Active	F05-12HG	Human	GST	Sf9 Cells	285-end (C491A)
FGFR2 (C491F), Active	F05-12IG	Human	GST	Sf9 Cells	285-end (C491F)
FGFR2 (C491S V564L), Active	F05-12MG	Human	GST	Sf9 Cells	285-end (C491S V564L)
FGFR2 (C491S), Active	F05-12JG	Human	GST	Sf9 Cells	285-end (C491S)
FGFR2 (E565A), Active	F05-12KG	Human	GST	Sf9 Cells	285-end (E565A)
FGFR2 (E565G), Active	F05-12CG	Human	GST	Sf9 Cells	285-end (E565G)
FGFR2 (K526E), Active	F05-12BG	Human	GST	Sf9 Cells	285-end (K526E)
FGFR2 (K641R), Active	F05-12DG	Human	GST	Sf9 Cells	285-end (K641R)
FGFR2 (K659N), Active	F05-12EG	Human	GST	Sf9 Cells	285-end (K659N)
FGFR2 (N549H), Active	F05-12G	Human	GST	Sf9 Cells	285-end (N549H)
FGFR2 (N549K), Active	F05-12NG	Human	GST	Sf9 Cells	285-end (N549K)
FGFR2 (R612T), Active	F05-12GG	Human	GST	Sf9 Cells	285-end (R612T)
FGFR2 (V564F), Active	F05-12FG	Human	GST	Sf9 Cells	285-end (V564F)
FGFR2 (V564I), Active	F05-12RG	Human	GST	Sf9 Cells	285-end (V564I)
FGFR2 (V564L), Active	F05-12LG	Human	GST	Sf9 Cells	285-end (V564L)
FGFR3 (G697C), Active	F06-12FG	Human	GST	Sf9 Cells	397-end (G697C)
FGFR3 (K650E), Active	F06-12CG	Human	GST	Sf9 Cells	397-end (K650E)
FGFR3 (K650M), Active	F06-12DG	Human	GST	Sf9 Cells	397-end (K650M)
FGFR3 (K650Q), Active	F06-12EG	Human	GST	Sf9 Cells	397-end (K650Q)
FGFR3 (V555M), Active	F06-12GG	Human	GST	Sf9 Cells	397-end (V555M)
FGFR4 (N535K), Active	F07-12G	Human	GST	Sf9 Cells	460-end (N535K)

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
FGFR4 (V550E), Active	F07-12BG	Human	GST	Sf9 Cells	460-end (V550E)
FGFR4 (V550L), Active	F07-12CG	Human	GST	Sf9 Cells	460-end (V550L)
FGFR4 (V550M), Active	F07-12DG	Human	GST	Sf9 Cells	460-end (V550M)
FIP1L1-PDGFR alpha, Active	P12-19BG	Human	GST	Sf9 Cells	1-265 (FIP1L1)
FLT3 (D835Y), Active	F12-12BG	Human	GST	Sf9 Cells	571-993 (D835Y)
FLT3 (F594_R595insR), Active	F12-12EG	Human	GST	Sf9 Cells	571-993 (F594_R595insR)
FLT3 (F594_R595insREY), Active	F12-12FG	Human	GST	Sf9 Cells	571-993 (F594_R595insREY)
FLT3 (F691L D835Y), Active	F12-12KG	Human	GST	Sf9 Cells	571-993 (F691L D835Y)
FLT3 (N676K), Active	F12-12IG	Human	GST	Sf9 Cells	571-993 (N676K)
FLT3 (R595_E596insEY), Active	F12-12GG	Human	GST	Sf9 Cells	571-993 (R595_E596insEY)
FLT3 (Y591-V592insVDFREYEYD D835Y), Active	F12-12JG	Human	GST	Sf9 Cells	571-993 (Y591-V592insVDFREY-EYD D835Y)
FLT3 (Y591-V592insVDFREYEYD F691L), Active	F12-12LG	Human	GST	Sf9 Cells	571-993 (Y591-V592insVDFREY-EYD F691L)
FLT3 (Y591-V592insVDFREYEYD), Active	F12-12HG	Human	GST	Sf9 Cells	571-993 (Y591-V592insVDFREY-EYD)
FLT3-ITD-NPOS, Active	F12-12DG	Human	GST	Sf9 Cells	571-993 (FLT3-ITD-NPOS)
FLT3-ITD-W51, Active	F12-12CG	Human	GST	Sf9 Cells	571-993 (FLT3-ITD-W51)
GOPC-ROS1, Active	R14-19BG	Human	GST	Sf9 Cells	1-419 (GOPC) - 1881-end (ROS1)
GSK3b (C199A), Active	G09-12H	Human	HIS	Sf9 Cells	Full-length (C199A)
HER2 (A775_G776insYVMA), Active	E27-13BG	Human	GST	Sf9 Cells	676-end (A775_G776insYVMA)
HER2 (d755-759), Active	E27-12BG	Human	GST	Sf9 Cells	676-end (d755-759)
HER2 (D769H), Active	E27-12GG	Human	GST	Sf9 Cells	676-end (D769H)
HER2 (D769Y), Active	E27-12HG	Human	GST	Sf9 Cells	676-end (D769Y)
HER2 (G776VC), Active	E27-13EG	Human	GST	Sf9 Cells	676-end (G776VC)
HER2 (I767M), Active	E27-12NG	Human	GST	Sf9 Cells	676-end (I767M)
HER2 (L755S), Active	E27-12EG	Human	GST	Sf9 Cells	676-end (L755S)
HER2 (L755W), Active	E27-12FG	Human	GST	Sf9 Cells	676-end (L755W)
HER2 (L785F), Active	E27-12LG	Human	GST	Sf9 Cells	676-end (L785F)
HER2 (P780_Y781insGSP), Active	E27-13DG	Human	GST	Sf9 Cells	676-end (P780_Y781insGSP)
HER2 (R896C), Active	E27-12KG	Human	GST	Sf9 Cells	676-end (R896C)
HER2 (T798I), Active	E27-12MG	Human	GST	Sf9 Cells	676-end (T798I)
HER2 (V777L), Active	E27-12IG	Human	GST	Sf9 Cells	676-end (V777L)
HER2 (V777_G778insCG), Active	E27-13CG	Human	GST	Sf9 Cells	676-end (V777_G778insCG)
HER2 (V842I), Active	E27-12JG	Human	GST	Sf9 Cells	676-end (V842I)
HER2 (V956R), Active	E27-12PH	Human	HIS	Sf9 Cells	677-1027 (V956R)
HER3 (E928G) Protein	E28-32AG	Human	GST	Sf9 Cells	684-1020 (E928G)
HER4 (R927Q), Active	E29-12G	Human	GST	Sf9 Cells	682-993 (R927Q)
InsR (R1201Q), Active	I08-122G	Human	GST	Sf9 Cells	980-1382 (R1201Q)
JAK1 (R724H), Active	J01-12BG	Human	GST	Sf9 Cells	438-end (R724H)
KDR (R1032Q) Protein	K01-12G	Human	GST	Sf9 Cells	789-end (R1032Q)
KIAA1549-BRAF (Kex15Bex9), Active	B08-19CG	Human	GST	Sf9 Cells	1320-1643 exon15(KIAA1549) - 381-end or exon9-18 (BRAF)

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
KIAA1549-BRAF (Kex16Bex9), Active	B08-19BG	Human	GST	Sf9 Cells	1320-1749 exon16 (KIAA1549) - 381-end or exon9-18 (BRAF)
KIF5B-ALK (Kex24Aex20), Active	A19-19MG	Human	GST	Sf9 Cells	1-921a.a., exons1-24 (KIF5B) - 1058a.a.-end, exons20-29 (ALK)
KIF5B-MET (Kex24Mex14), Active	M52-19BG	Human	GST	Sf9 Cells	1-921a.a., exons 1-24 (KIF5B) 963 a.a. - end, exons 14-21 (MET)
KIF5B-RET (Kex15Rex14), Active	R02-19FG	Human	GST	Sf9 Cells	1- 575 a.a., exons1-15 (KIF5B) – 713 a.a. – end, exons 14-21 (RET)
KIF5B-RET (V804L), Active	R02-19HG	Human	GST	Sf9 Cells	1- 575 a.a., exons1-15 (KIF5B) – 713 a.a. – end, exons 14-21 (RET (V804L))
KIF5B-RET (V804M), Active	R02-19GG	Human	GST	Sf9 Cells	1- 575 a.a., exons1-15 (KIF5B) - 713 a.a. – end, exons 14-21 (RET (V804M))
KLC1-ALK (Kex8Aex20), Active	A19-19IG	Human	GST	Sf9 Cells	1-420a.a., exons1-8 (KLC1)-1058a.a.-end, exons20-29 (ALK)
LRRK2 (D1994A), Active	L10-12FG	Human	GST	Sf9 Cells	968-end (D1994A)
LRRK2 (G2019S), Active	L10-12GG	Human	GST	Sf9 Cells	968-end (G2019S)
LRRK2 (G2385R), Active	L10-12IG	Human	GST	Sf9 Cells	968-end (G2385R)
LRRK2 (I1371V), Active	L10-12JG	Human	GST	Sf9 Cells	968-end (I1371V)
LRRK2 (I2020T), Active	L10-12HG	Human	GST	Sf9 Cells	968-end (I2020T)
LRRK2 (R1441C), Active	L10-12G	Human	GST	Sf9 Cells	968-end (R1441C)
LRRK2 (R1441G), Active	L10-12BG	Human	GST	Sf9 Cells	968-end (R1441G)
LRRK2 (R1441H), Active	L10-12CG	Human	GST	Sf9 Cells	968-end (R1441H)
LRRK2 (Y1699C), Active	L10-12DG	Human	GST	Sf9 Cells	968-end (Y1699C)
LRRK2 (Y1699G), Active	L10-12EG	Human	GST	Sf9 Cells	968-end (Y1699G)
MELK (T460M), Active	M50-12G	Human	GST	Sf9 Cells	1-550 (T460M)
MET (D1213H), Active	M52-12F2G	Human	GST	Sf9 Cells	956-end (D1213H)
MET (D1228A), Active	M52-12ZG	Human	GST	Sf9 Cells	956-end (D1228A)
MET (D1228G), Active	M52-12B2G	Human	GST	Sf9 Cells	956-end (D1228G)
MET (D1228H), Active	M52-12HG	Human	GST	Sf9 Cells	956-end (D1228H)
MET (D1228N), Active	M52-12IG	Human	GST	Sf9 Cells	956-end (D1228N)
MET (D1228V), Active	M52-12C2G	Human	GST	Sf9 Cells	956-end (D1228V)
MET (D1228Y), Active	M52-12D2G	Human	GST	Sf9 Cells	956-end (D1228Y)
MET (Del Ex14), Active	M52-12PG	Human	GST	Sf9 Cells	956-end (Del Ex14)
MET (F1200I), Active	M52-12GG	Human	GST	Sf9 Cells	956-end (F1200I)
MET (G1090A), Active	M52-12E2G	Human	GST	Sf9 Cells	956-end (G1090A)
MET (G1163R), Active	M52-12XG	Human	GST	Sf9 Cells	956-end (G1163R)
MET (H1094L), Active	M52-12VG	Human	GST	Sf9 Cells	956-end (H1094L)
MET (H1094Y), Active	M52-12WG	Human	GST	Sf9 Cells	956-end (H1094Y)
MET (K1244R), Active	M52-12QG	Human	GST	Sf9 Cells	956-end (K1244R)
MET (L1195F), Active	M52-12YG	Human	GST	Sf9 Cells	956-end (L1195F)
MET (L1195V), Active	M52-12TG	Human	GST	Sf9 Cells	956-end (L1195V)
MET (M1250I), Active	M52-12SG	Human	GST	Sf9 Cells	956-end (M1250I)
MET (M1250T), Active	M52-12BG	Human	GST	Sf9 Cells	956-end (M1250T)

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
<b>MET (P991S), Active</b>	M52-12CG	Human	GST	Sf9 Cells	956-end (P991S)
<b>MET (R970C), Active</b>	M52-12OG	Human	GST	Sf9 Cells	956-end (R970C)
<b>MET (R1227K), Active</b>	M52-12G2G	Human	GST	Sf9 Cells	956-end (R1227K)
<b>MET (T1173I), Active</b>	M52-12FG	Human	GST	Sf9 Cells	956-end (T1173I)
<b>MET (T992I), Active</b>	M52-12DG	Human	GST	Sf9 Cells	956-end (T992I)
<b>MET (V1092I), Active</b>	M52-12EG	Human	GST	Sf9 Cells	956-end (V1092I)
<b>MET (Y1230A), Active</b>	M52-12JG	Human	GST	Sf9 Cells	956-end (Y1230A)
<b>MET (Y1230C), Active</b>	M52-12KG	Human	GST	Sf9 Cells	956-end (Y1230C)
<b>MET (Y1230D), Active</b>	M52-12LG	Human	GST	Sf9 Cells	956-end (Y1230D)
<b>MET (Y1230H), Active</b>	M52-12MG	Human	GST	Sf9 Cells	956-end (Y1230H)
<b>MET (Y1230S), Active</b>	M52-12UG	Human	GST	Sf9 Cells	956-end (Y1230S)
<b>MET (Y1235D), Active</b>	M52-12NG	Human	GST	Sf9 Cells	956-end (Y1235D)
<b>NPM1-ALK (F1174L), Active</b>	A19-19EG	Human	GST	Sf9 Cells	1-117 (NPM1) 1058-end (F1174L) (ALK)
<b>NPM1-ALK, Active</b>	A19-19BG	Human	GST	Sf9 Cells	1-117 (NPM1) - 1058-end (ALK)
<b>p70S6Kb (C150T), Active</b>	R22-12G	Human	GST	Sf9 Cells	Full length (C150T)
<b>PAK2 (Y443N), Active</b>	P03-12G	Human	GST	Sf9 Cells	Full length (Y443N)
<b>PDGFR alpha (D842V), Active</b>	P12-12BG	Human	GST	Sf9 Cells	550-end (D842V)
<b>PDGFR alpha (E675G Y676C), Active</b>	P12-12EG	Human	GST	Sf9 Cells	550-end (E675G Y676C)
<b>PDGFR alpha (G680R), Active</b>	P12-12DG	Human	GST	Sf9 Cells	550-end (G680R)
<b>PDGFR alpha (T674I), Active</b>	P12-12CG	Human	GST	Sf9 Cells	550-end (T674I)
<b>PFKFB3 (S269A), Active</b>	P323-10BG	Human	GST	Sf9 Cells	Full length (S269A)
<b>PI3K (p110 alpha (E542K)/p85 alpha), Active</b>	P27-122DH	Human	HIS	Sf9 Cells	Full Length (p110 alpha (E542K)/p85 alpha)
<b>PI3K (p110 alpha (E545K)/p85 alpha), Active</b>	P27-152H	Human	HIS	Sf9 Cells	Full Length (p110 alpha (E545K)/p85 alpha)
<b>PI3K (p110 alpha (E545K)/p85 alpha), Active</b>	P27-152G	Human	GST	Sf9 Cells	Full length (p110 alpha (E545K)/p85 alpha)
<b>PI3K (p110 alpha (N345K)/p85 alpha), Active</b>	P27-122CH	Human	HIS	Sf9 Cells	Full Length (p110 alpha (N345K)/p85 alpha)
<b>PI3K (p110 alpha(H1047R)/p85 alpha), Active</b>	P27-122EH	Human	HIS	Sf9 Cells	Full length (p110 alpha(H1047R)/p85 alpha)
<b>PI3K (p110 delta (E1021K)/p85 alpha), Active</b>	P30-122H	Human	HIS	Sf9 Cells	Full-Length (p110 delta (E1021K)/p85 alpha)
<b>PKD2 (PRKD2) (G848E), Active</b>	P76-12BG	Human	GST	Sf9 Cells	Full Length (G848E)
<b>PKD2 (PRKD2) (G870E), Active</b>	P76-12CG	Human	GST	Sf9 Cells	Full Length (G870E)
<b>PLK4 (G95L), Active</b>	P44-12G	Human	GST	Sf9 Cells	1-836 (G95L)
<b>PRKACA (L206R), Active</b>	P51-19CG	Human	GST	Sf9 Cells	1-351 (L206R)
<b>PTC1 (CCDC6-RET), Active</b>	R02-19BG	Human	GST	Sf9 Cells	CCDC6 (1-101)-RET (713-end)
<b>PTC2 (PRKAR1A-RET), Active</b>	R02-19CG	Human	GST	Sf9 Cells	PRKAR1A (1-236)-RET (713-end)
<b>PTC3 (NCOA4-RET), Active</b>	R02-19DG	Human	GST	Sf9 Cells	NCOA4 (1-238)-RET (713-end)
<b>RAF1 (EE), Active</b>	R01-13G	Human	GST	Sf9 Cells	Y340E-Y341E, 306-end
<b>RAF1 (R391W), Active</b>	R01-12BG	Human	GST	Sf9 Cells	307-end (R391W)

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
RET (A883F), Active	R02-12MG	Human	GST	Sf9 Cells	658-end (A883F)
RET (E762Q), Active	R02-12FG	Human	GST	Sf9 Cells	658-end (E762Q)
RET (G691S), Active	R02-12DG	Human	GST	Sf9 Cells	658-end (G691S)
RET (G810R), Active	R02-12VG	Human	GST	Sf9 Cells	658-end (G810R)
RET (G810S M918T), Active	R02-12WG	Human	GST	Sf9 Cells	658-end (G810S M918T)
RET (G810V), Active	R02-12QG	Human	GST	Sf9 Cells	658-end (G810V)
RET (L790F), Active	R02-12LG	Human	GST	Sf9 Cells	658-end (L790F)
RET (M918T), Active	R02-12JG	Human	GST	Sf9 Cells	658-end (M918T)
RET (M918T), Active	R02-12JH	Human	HIS	Sf9 Cells	658-end (M918T)
RET (R749T), Active	R02-12EG	Human	GST	Sf9 Cells	658-end (R749T)
RET (R813Q), Active	R02-12HG	Human	GST	Sf9 Cells	658-end (R813Q)
RET (R912P), Active	R02-12PG	Human	GST	Sf9 Cells	658-end (R912P)
RET (S891A), Active	R02-12IG	Human	GST	Sf9 Cells	658-end (S891A)
RET (S904A), Active	R02-12OG	Human	GST	Sf9 Cells	658-end (S904A)
RET (S904A), Active	R02-12OH	Human	HIS	Sf9 Cells	658-end (S904A)
RET (S904F), Active	R02-12NG	Human	GST	Sf9 Cells	658-end (S904F)
RET (S904F), Active	R02-12NH	Human	HIS	Sf9 Cells	658-end (S904F)
RET (V738A), Active	R02-12RG	Human	GST	Sf9 Cells	658-end (V738A)
RET (V778I), Active	R02-12KG	Human	GST	Sf9 Cells	658-end (V778I)
RET (V804L), Active	R02-12BG	Human	GST	Sf9 Cells	658-end (V804L)
RET (V804M G810S), Active	R02-12UG	Human	GST	Sf9 Cells	658-end (V804M G810S)
RET (V804M), Active	R02-12GG	Human	GST	Sf9 Cells	658-end (V804M)
RET (Y791F), Active	R02-12CG	Human	GST	Sf9 Cells	658-end (Y791F)
RET (Y806C), Active	R02-12SG	Human	GST	Sf9 Cells	658-end (Y806C)
RET (Y806N), Active	R02-12TG	Human	GST	Sf9 Cells	658-end (Y806N)
RNF130-BRAF (Rex3Bex9), Active	B08-19GG	Human	GST	Sf9 Cells	1-231 exon3 (RNF130) - 381-end or exon 9-18 (BRAF)
ROS1 (G2032R), Active	R14-12BG	Human	GST	Sf9 Cells	1881-end (G2032R)
ROS1 (G2101A), Active	R14-12EG	Human	GST	Sf9 Cells	1881-end (G2101A)
ROS1 (G2101C), Active	R14-12DG	Human	GST	Sf9 Cells	1881-end (G2101C)
ROS1(L2086F), Active	R14-12CG	Human	GST	Sf9 Cells	1881-end (L2086F)
RSK2 (I416V), Active	R17-12BG	Human	GST	Sf9 Cells	Full Length (I416V)
RSK2 (L608F), Active	R17-12CG	Human	GST	Sf9 Cells	Full Length (L608F)
RSK2 (Y483C), Active	R17-12DG	Human	GST	Sf9 Cells	Full Length (Y483C)
SRGAP3-BRAF (Sex12Bex9), Active	B08-19FG	Human	GST	Sf9 Cells	1-513 exon12 (SRGAP3) 381-end or exon 9-18 (BRAF)
STRN-ALK (Sex3Aex20), Active	A19-19JG	Human	GST	Sf9 Cells	1-137 (STRN)-1058-end (ALK)
TECR-PKN1(Tex1Pex10), Active	P70-19BG	Human	GST	Sf9 Cells	1-5a.a., exons1 (TECR) - 473a.a.-end, exons10-22 (PKN1)
TFG-ALK (Tex4Aex20), Active	A19-19KG	Human	GST	Sf9 Cells	1-138a.a., exons1-4 (TFG) - 1058a.a.-end, exons20-29 (ALK)
TFG-ALK (Tex5Aex20), Active	A19-19CG	Human	GST	Sf9 Cells	1-193 (TFG)-1058-end (ALK)
TFG-MET (Tex5Mex15), Active	M52-19G	Human	GST	Sf9 Cells	1-193a.a., exons1-5 (TFG) - 1010a.a.-end, exons15-21 (MET)

# Mutant Kinases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
TFG-TRKA (TRK-T3), Active	N16-19CG	Human	GST	Sf9 Cells	1-193 (TFG)-399-end (TRKA)
TGFBR1 (T204D), Active	T07-13G	Human	GST	Sf9 Cells	80-end (T204D)
TIE2 (A1124V), Active	T04-12GG	Human	GST	Sf9 Cells	771-end (A1124V)
TIE2 (L914F), Active	T04-12IG	Human	GST	Sf9 Cells	771-end (L914F)
TIE2 (P883A), Active	T04-12CG	Human	GST	Sf9 Cells	771-end (P883A)
TIE2 (R849W), Active	T04-12BG	Human	GST	Sf9 Cells	771-end (R849W)
TIE2 (R915C), Active	T04-12JG	Human	GST	Sf9 Cells	771-end (R915C)
TIE2 (Y1108F), Active	T04-12FG	Human	GST	Sf9 Cells	771-end (Y1108F)
TIE2 (Y897C), Active	T04-12DG	Human	GST	Sf9 Cells	771-end (Y897C)
TIE2 (Y897H R915C), Active	T04-12KG	Human	GST	Sf9 Cells	771-end (Y897H R915C)
TIE2 (Y897H), Active	T04-12HG	Human	GST	Sf9 Cells	771-end (Y897H)
TIE2 (Y897S), Active	T04-12EG	Human	GST	Sf9 Cells	771-end (Y897S)
TPM1-ALK, Active	A19-19LG	Human	GST	Sf9 Cells	TPM1 (1-257)-ALK (1058-end)
TPM3-ALK, Active	A19-19DG	Human	GST	Sf9 Cells	1-258 (TPM3) - 1058-end (ALK)
TPM3-PDGFR beta, Active	P13-19DG	Human	GST	Sf9 Cells	1-258 (TPM3) - 528-end (PDGFRB)
TPM3-ROS1, Active	R14-19CG	Human	GST	Sf9 Cells	1-258 (TPM3) - 1881-end (ROS1)
TPM3-TRKA, Active	N16-19DG	Human	GST	Sf9 Cells	1-258 (TPM3) - 399-end (TRKA)
TPR-TRKA (TRK-T1), Active	N16-19BG	Human	GST	Sf9 Cells	1-191(TPR) - 399-end (TRKA)
TRKA (A608D), Active	N16-12DG	Human	GST	Sf9 Cells	440-end (A608D)
TRKA (F589L), Active	N16-12EG	Human	GST	Sf9 Cells	440-end (F589L)
TRKA (G595R A608D), Active	N16-12FG	Human	GST	Sf9 Cells	440-end (G595R A608D)
TRKA (G595R G667A), Active	N16-12JG	Human	GST	Sf9 Cells	440-end (G595R G667A)
TRKA (G595R G667C), Active	N16-12KG	Human	GST	Sf9 Cells	440-end (G595R G667C)
TRKA (G595R G667S), Active	N16-12LG	Human	GST	Sf9 Cells	440-end (G595R G667S)
TRKA (G595R L657M), Active	N16-12GG	Human	GST	Sf9 Cells	440-end (G595R L657M)
TRKA (G595R), Active	N16-12BG	Human	GST	Sf9 Cells	440-end (G595R)
TRKA (G667A), Active	N16-12MG	Human	GST	Sf9 Cells	440-end (G667A)
TRKA (G667C), Active	N16-12CG	Human	GST	Sf9 Cells	440-end (G667C)
TRKA (G667S), Active	N16-12NG	Human	GST	Sf9 Cells	440-end (G667S)
TRKA (L657M), Active	N16-12HG	Human	GST	Sf9 Cells	440-end (L657M)
TRKC (G623E), Active	N18-12BH	Human	HIS	Sf9 Cells	507-end (G623E)
TRKC (G623R L686M), Active	N18-12EH	Human	HIS	Sf9 Cells	507-end (G623R L686M)
TRKC (G623R), Active	N18-12CH	Human	HIS	Sf9 Cells	507-end (G623R)
TRKC (G696A), Active	N18-12FH	Human	HIS	Sf9 Cells	454-end (G696A)
TRKC (G696C), Active	N18-12GH	Human	HIS	Sf9 Cells	454-end (G696C)
TRKC (L686M), Active	N18-12DH	Human	HIS	Sf9 Cells	507-end (L686M)
YES1 (T348I), Active	Y01-12BG	Human	GST	Sf9 Cells	Full Length (T348I)
ZAP70 (Y319F), Active	Z02-12BG	Human	GST	Sf9 Cells	Full Length (Y319F)

# Phosphatases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
CaMKPase Protein	P06-34G	Human	GST	E.coli	Full Length
CDC25A, Active	C04-20G	Human	GST	Sf9 Cells	Full Length
CDC25B, Active	C04-20BG	Human	GST	Sf9 Cells	Full Length
CDC25C, Active	C04-20CG	Human	GST	Sf9 Cells	Full Length
DUSP22 (MKPX), Active	D22-20G	Human	GST	E.coli	Full Length
Lambda Protein Phosphatase	L01-20U	Bacte- riophage Lambda		E.coli	Full length
PHLPP2 Protein	P72-21G	Human	GST	E.coli	766-1043
PP1A, Active	P13-20G	Human	GST	Sf9 Cells	Full Length
PP1A, Active	P13-20GH	Human	GST	E.coli	Full length
PP1B, Active	P14-20G	Human	GST	Sf9 Cells	Full Length
PP2A alpha/PPP2R1A Complex, Active	P16-20BH	Human	HIS	Sf9 Cells	Full Length
PP2Aα/PPP2R1A Complex, Active	P16-20BG	Human	GST	Sf9 Cells	Full length
PP2Abeta Protein	P17-34G	Human	GST	Sf9 Cells	Full Length
PP2Calpha, Active	P02-20G	Human	GST	Sf9 Cells	Full Length
PP2Calpha, Active	P02-20BG	Human	GST	E.coli	Full Length
PP2Cgamma, Active	P07-20G	Human	GST	Sf9 Cells	Full Length
PTEN Protein	P23-34G	Human	GST	E.coli	Full Length
PTEN, Active	P23-20G	Human	GST	E.coli	Full Length
PTEN, Active	T03-20H	Human	HIS	Sf9 Cells	Full length
PTPN1 (PTP1B), Active	P28-20G	Human	GST	Sf9 Cells	Full Length
PTPN11 (SHP2), Active	P38-21G	Human	GST	E.coli	246-593
PTPN11 (SHP2), Active	P38-20G	Human	GST	Sf9 Cells	Full Length
PTPN11 (T253M Q257L), Active	P38-22G	Human	GST	Sf9 Cells	Full Length
PTPN12 (PTP-PEST), Active	P39-21G	Human	GST	E.coli	1-355
PTPN13 (FAP-1), Active	P40-21G	Human	GST	E.coli	2169-2485
PTPN2 (TC-PTP), Active	P29-20G	Human	GST	E.coli	Full Length
PTPN6 (SHP1), Active	P33-20G	Human	GST	E.coli	Full Length
PTPN6 (SHP1), Active	P33-20H	Human	HIS	E.coli	Full Length
PTPN7 (LC-PTP), Active	P34-20G	Human	GST	E.coli	Full Length
PTPRA (LRP), Active	P48-21G	Human	GST	E.coli	174-802
PTPRB, Active	P49-21G	Human	GST	Sf9 Cells	1552-end
PTPRC (CD45), Active	P50-21H	Human	HIS	Sf9 Cells	604-end
PTPRC (CD45), Active	P50-21BG	Mouse	GST	E.coli	592-end
PTPRE, Active	P52-21G	Human	GST	Sf9 Cells	75-end
PTPRF (LAR), Active	P53-21G	Human	GST	E.coli	1275-1897
PTPRJ (CD148), Active	P56-21G	Human	GST	Sf9 Cells	Full Length
PTPRM (PTP- mu) Protein	P58-35G	Human	GST	Sf9 Cells	764-1452
PTPRS (PTP- sigma), Active	P64-21G	Human	GST	E.coli	883-1210
SHIP1, Active	P08-20G	Human	GST	Sf9 Cells	Full Length
SHIP2, Active	P09-20G	Human	GST	Sf9 Cells	Full Length
WIP1, Active	P04-20G	Human	GST	Sf9 Cells	Full length

# Epigenetic Enzymes

## HISTONE Deacetylases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
HDAC1, Active	H83-30G	Human	GST	Sf9 Cells	Full Length
HDAC10, Active	H92-31G	Human	GST	Sf9 Cells	1-482
HDAC11, Active	H93-30G	Human	GST	Sf9 Cells	Full Length
HDAC11, Active	H93D-30G	Dog	GST	Sf9 Cells	Full length
HDAC11, Active	H93C-30G	Macaca fascicularis	GST	Sf9 Cells	Full length
HDAC11, Active	H93M-30G	Mouse	GST	Sf9 Cells	Full length
HDAC11, Active	H93A-30G	Rat	GST	Sf9 Cells	Full length
HDAC2, Active	H84-30G	Human	GST	Sf9 Cells	Full Length
HDAC3, Active	H85-30G	Human	GST	Sf9 Cells	Full Length
HDAC3/NCOR2, Active	H85-38G	Human	GST	Sf9 Cells	HDAC3 (full-Length)/ NCOR2 (237-489)
HDAC4, Active	H86-31G	Human	GST	Sf9 Cells	612-end
HDAC4, Active	H86-31BG	Human	GST	Sf9 Cells	101-end
HDAC5, Active	H87-31G	Mouse	GST	Sf9 Cells	617-end
HDAC6, Active	H88-30G	Human	GST	Sf9 Cells	Full Length
HDAC6, Active	H88M-30G	Mouse	GST	Sf9 Cells	Full length
HDAC6, Active	H88A-30G	Rat	GST	Sf9 Cells	Full length
HDAC6, Active	H88D-30G	Dog	GST	Sf9 Cells	Full length
HDAC6, Active	H88C-30G	Macaca fascicularis	GST	Sf9 Cells	Full length
HDAC7, Active	H89-31G	Human	GST	Sf9 Cells	501-end
HDAC8, Active	H90-30H	Human	HIS	Sf9 Cells	Full Length
HDAC8, Active	H90-30G	Human	GST	Sf9 Cells	Full Length
HDAC9, Active	H91-31G	Human	GST	Sf9 Cells	548-end
SIRT1, Active	S35-31H	Human	HIS	Sf9 Cells	193-end
SIRT1, Active	S35-31G	Human	GST	Sf9 Cells	193-end
SIRT2, Active	S36-30H	Human	HIS	Sf9 Cells	Full Length
SIRT3, Active	S37-38G	Human	GST	Sf9 Cells	47-end
SIRT5, Active	S39-30H	Human	HIS	Sf9 Cells	Full Length
SIRT6, Active	S40-31H	Human	HIS	Sf9 Cells	23-end
SIRT7, Active	S41-30H	Human	HIS	Sf9 Cells	Full Length

# Acetyl and Methyltransferases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
ASH2L Protein	A372-30BG	Human	GST	E.coli	Full Length
COMT, Active	C339-381G	Human	GST	Sf9 Cells	27-271
DNMT1 Protein	D351-31G	Mouse	GST	Sf9 Cells	766-end
DNMT3A, Active	D353-380G	Human	GST	Sf9 Cells	Full Length
DNMT3B, Active	D353-380BG	Human	GST	Sf9 Cells	Full Length
DNMT3L, Active	D353-380CG	Human	GST	Sf9 Cells	Full Length
DOT1L (KMT4), Active	D344-381G	Human	GST	Sf9 Cells	1-435
DPY30 Protein	D369-30H	Human	HIS	E.coli	Full Length
EED Protein	E310-30H	Human	HIS	E.coli	Full Length
EHMT1 (KMT1D), Active	E317-381G	Human	GST	Sf9 Cells	895-end
EHMT2, Active	E318-381G	Human	GST	Sf9 Cells	807-end
EZH1 Protein	E395-30G	Human	GST	Sf9 Cells	Full Length
EZH2 (Y641C) Protein	E396-32BG	Human	GST	Sf9 Cells	Full Length
EZH2 (Y641H) Protein	E396-32DG	Human	GST	Sf9 Cells	Full Length
EZH2 (Y641N) Protein	E396-32EG	Human	GST	Sf9 Cells	Full Length
EZH2 (Y641S) Protein	E396-32FG	Human	GST	Sf9 Cells	Full Length
EZH2 Protein	E396-30G	Human	GST	Sf9 Cells	Full Length
HEMK1 Protein	H311-30G	Human	GST	Sf9 Cells	Full Length
KAT2A (GCN5), Active	K311-381G	Human	GST	Sf9 Cells	323-end
KAT2B (PCAF), Active	K311-381BG	Human	GST	Sf9 Cells	431-end
KAT3A (CREBBP) Protein	K312-31G	Human	GST	Sf9 Cells	518-1207
KAT3A (CREBBP), Active	K312-381G	Human	GST	Sf9 Cells	1319-1710
KAT3B (EP300) Protein	K312-31BG	Human	GST	Sf9 Cells	532-1153; contains the catalytic domain
KAT5 (TIP60), Active	K314-380G	Human	GST	Sf9 Cells	Full Length
KAT6B (MYST4), Active	K315-381BG	Human	GST	Sf9 Cells	431-end
KAT7 (MYST2), Active	K316-380G	Human	GST	Sf9 Cells	Full Length
KAT8 (MYST1) Protein	K317-31G	Human	GST	Sf9 Cells	2-467
KAT9 (ELP3) Protein	K318-30G	Human	GST	Sf9 Cells	Full Length
LCMT1, Active	L321-380G	Human	GST	Sf9 Cells	Full Length
LCMT2 Protein	L322-30G	Human	GST	Sf9 Cells	Full Length
MEP50 (WDR77) Protein	W329-30G	Human	GST	Sf9 Cells	Full Length
METTL1 Protein	M321-30G	Human	GST	Sf9 Cells	Full Length
METTL11A Protein	M331-30G	Human	GST	Sf9 Cells	Full Length
METTL12, Active	M332-30G	Human	GST	Sf9 Cells	Full length
METTL14 Protein	M334-30G	Human	GST	Sf9 Cells	Full Length
METTL16, Active	M336-380G	Human	GST	Sf9 Cells	Full length
METTL2A Protein	M322-30G	Human	GST	Sf9 Cells	Full Length
METTL3 Protein	M323-30G	Human	GST	Sf9 Cells	Full Length
METTL3/METTL14, Active	M323-380G	Human	GST	Sf9 Cells	Full Length
MGMT Protein	M354-30H	Human	HIS	Sf9 Cells	Full Length
MLL1 (KMT2A), Active	M341-381G	Human	GST	Sf9 Cells	3738-end
MLL2 (KMT2D), Active	M342-381G	Human	GST	Sf9 Cells	5317-end
MLL3 (KMT2C), Active	M343-381G	Human	GST	Sf9 Cells	4689-end
MLL5 (KMT2E), Active	M345-381G	Human	GST	Sf9 Cells	1-548

# Acetyl and Methyltransferases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
N6AMT2, Active	N355-380BH	Human	HIS	Sf9 Cells	Full Length
NNMT (Dog), Active	N330D-380G	Canis lupus	GST	Sf9 Cells	Full Length
NNMT, Active	N330-380G	Human	GST	Sf9 Cells	Full Length
PCAF Protein	P09-31G	Human	GST	E.coli	431-end
PCMT1 Protein	P314-30G	Human	GST	Sf9 Cells	2-end
PNMT, Active	P331-380G	Human	GST	Sf9 Cells	Full Length
PRDM2 (KMT8) Protein	P333-31G	Human	GST	Sf9 Cells	1-340
PRMT1, Active	P365-380G	Human	GST	Sf9 Cells	Full Length
PRMT2 Protein	P365-30BG	Human	GST	Sf9 Cells	Full Length
PRMT3, Active	P365-380CG	Human	GST	Sf9 Cells	Full Length
PRMT4, Active	P365-380DG	Human	GST	Sf9 Cells	Full Length
PRMT5 Protein	P365-30EG	Human	GST	Sf9 Cells	Full Length
PRMT6, Active	P365-380FG	Human	GST	Sf9 Cells	Full Length
PRMT7 Protein	P365-30GG	Human	GST	Sf9 Cells	Full Length
PRMT8, Active	P365-380HG	Human	GST	Sf9 Cells	Full Length
RBBP5 Protein	R315-30H	Human	HIS	E.coli	Full Length
SETD1A (KMT2F), Active	S342-381G	Human	GST	Sf9 Cells	1418-end
SETD2 (KMT3A), Active	S343-381G	Human	GST	Sf9 Cells	1425-1717
SETD4 Protein	S353-30G	Human	GST	Sf9 Cells	Full Length
SETD6, Active	S355-380G	Human	GST	Sf9 Cells	Full Length
SETD7, Active	S344-380G	Human	GST	Sf9 Cells	Full Length
SETD8 Protein	S345-30G	Human	GST	Sf9 Cells	Full Length
SETDB1 (KMT1E) Protein	S346-31G	Human	GST	Sf9 Cells	715-end
SETDB2 Protein	S346-30BG	Human	GST	Sf9 Cells	Full Length
SETMAR Protein	S347-30G	Human	GST	Sf9 Cells	Full Length
SMYD2 (KMT3C), Active	S348-380BG	Human	GST	Sf9 Cells	Full Length
SMYD3, Active	S348-380CG	Human	GST	Sf9 Cells	Full Length
SUV39H1 (KMT1A), Active	S350-380G	Human	GST	Sf9 Cells	Full Length
SUV39H2 (KMT1B), Active	S350-380BG	Human	GST	Sf9 Cells	Full Length
SUV420H1 (KMT5B), Active	S351-380G	Human	GST	Sf9 Cells	Full Length
TRDMT1 (DNMT2) Protein	T352-30G	Human	GST	Sf9 Cells	Full Length
WDR5 Protein	W325-30H	Human	HIS	E.coli	Full Length
WHSC1 (NSD2), Active	N377-381G	Human	GST	Sf9 Cells	911-end
WHSC1L1 (NSD3) Protein	N378-31G	Human	GST	Sf9 Cells	1038-end
WTAP Protein	W369-30G	Human	GST	Sf9 Cells	Full Length

# Deacetylase & Demethylase Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
JMJD6 Protein	J453-30G	Human	GST	Sf9 Cells	Full Length
KDM1A Protein	K421-31G	Human	GST	Sf9 Cells	172-end
KDM1B Protein	K421-30BG	Mouse	GST	Sf9 Cells	Full Length
KDM2A Protein	K422-31G	Human	GST	Sf9 Cells	1-748

# Deacetylase & Demethylase Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
KDM2B Protein	K422-30BG	Human	GST	Sf9 Cells	Full Length
KDM4A Protein	K424-31G	Human	GST	Sf9 Cells	1-886
KDM4B Protein	K424-31BG	Mouse	GST	Sf9 Cells	1-725
KDM4C Protein	K424-31CG	Human	GST	Sf9 Cells	1-460
KDM5B Protein	K425-31BG	Mouse	GST	Sf9 Cells	1-753
KDM5C Protein	K425-31CG	Human	GST	Sf9 Cells	1-671
PPME1 Protein	P332-30G	Human	GST	Sf9 Cells	2-end
SIRT1 Protein	S35-31EG	Human	GST	E.coli	193-end
SIRT2 Protein	S36-30EG	Human	GST	E.coli	Full Length
SIRT3 Protein	S37-30EG	Human	GST	E.coli	Full Length
SIRT4 Protein	S38-30EG	Human	GST	E.coli	Full Length
SIRT5 Protein	S39-30EG	Human	GST	E.coli	Full Length
SIRT6 Protein	S40-30EG	Human	GST	E.coli	Full Length

## Bromodomains

Product Name	Cat# No	Species	Tag	Expression System	Sequence
BRD2 (BD2), Active	B16-31H	Human	HIS	E. Coli	339-459
BRD4 (BD1), Active	B18-31H	Human	HIS	E. Coli	49-170
BRD4 (BD2) , Active	B18-38H	Human	HIS	E. Coli	342-460

# Ubiquitin

## Ubiquitin Enzymes - E1

Product Name	Cat# No	Species	Tag	Expression System	Sequence
SAE1/UBA2 Complex, Active	U208-380G	Human	His/GST	Sf9 Cells	Full Length
UBA1 (UBE1), Active	U201-380G	Human	GST	Sf9 Cells	Full Length
UBA6, Active	U206-380G	Human	GST	Sf9 Cells	Full Length
UBA7(UBE1L), Active	U207-380G	Human	GST	Sf9 Cells	Full Length
UBA2 (SAE2) Protein	U202-30G	Human	GST	Sf9 Cells	Full Length
UBA3 (UBE1C) Protein	U203-30G	Human	GST	Sf9 Cells	Full Length
UBLE1A (SAE1) Protein	U208-30G	Human	GST	Sf9 Cells	Full Length
NAE1/UBA3 Complex, Active	U203-380G	Human	His/GST	Sf9 Cells	2-end

# Ubiquitin Enzymes - E2

Product Name	Cat# No	Species	Tag	Expression System	Sequence
UBE2A, Active	U210-380H	Human	HIS	E.coli	Full Length
UBE2C, Active	U212-380H	Human	HIS	E.coli	Full Length
UBE2D1 (UBCH5A), Active	U213-380H	Human	HIS	E.coli	Full Length
UBE2D2 (UBC4), Active	U214-380H	Xenopus	HIS	E.coli	Full Length
UBE2D3 (UBCH5C), Active	U215-380H	Human	HIS	E.coli	Full Length
UBE2E1 (UBCH6), Active	U217-380H	Human	HIS	E.coli	2-end
UBE2E3 (UBCH9), Active	U219-380H	Human	HIS	E.coli	Full Length
UBE2G1 (UBC7), Active	U221-380H	Human	HIS	E.coli	Full Length
UBE2I (UBC9), Active	U224-380H	Human	HIS	E.coli	Full Length
UBE2K (UBC1), Active	U227-380H	Human	HIS	E.coli	Full Length
UBE2L6 (UBCH8), Active	U229-380H	Human	HIS	E.coli	Full Length
UBE2N (UBC13), Active	U231-380H	Human	HIS	E.coli	2-end
UBE2O, Active	U232-380G	Human	GST	Sf9 Cells	Full Length
UBE2B Protein	U211-30H	Human	HIS	E.coli	Full Length
UBE2F Protein	U220-30H	Human	HIS	E.coli	Full Length
UBE2H Protein	U223-30H	Human	HIS	E.coli	Full Length
UBE2J1 Protein	U225-30G	Human	GST	Sf9 Cells	Full Length
UBE2L3 (UBCH7) Protein	U228-30H	Human	HIS	E.coli	Full Length
UBE2R1 Protein	U234-30H	Human	HIS	E.coli	Full Length
UBE2R2 (UBC3B) Protein	U235-30H	Human	HIS	E.coli	2-end
UBE2U Protein	U237-30G	Human	GST	Sf9 Cells	Full Length
UBE2V2 Protein	U239-30H	Human	HIS	E.coli	2-end
UBE2W (UBC16) Protein	U240-30G	Human	GST	E.coli	Full Length
UBE2Z (USE1) Protein	U241-31H	Human	HIS	E.coli	109-end

# Ubiquitin Enzymes - E3

Product Name	Cat# No	Species	Tag	Expression System	Sequence
BIRC3, Active	B280-380G	Human	GST	Sf9 Cells	Full Length
BIRC7, Active	B281-380G	Human	GST	Sf9 Cells	Full Length
HERC4, Active	H265-381G	Human	GST	Sf9 Cells	642-end
MGRN1, Active	M287-380G	Human	GST	Sf9 Cells	2-end
RanBP2, Active	R230-381H	Human		E.coli	2553-2838
RNF34 (CARP), Active	R296-380G	Human	GST	Sf9 Cells	Full Length
RNF34L (CARP2), Active	R297-380G	Human	GST	Sf9 Cells	Full Length
TRIM37, Active	T292-380G	Human	GST	Sf9 Cells	Full Length
WWP2, Active	W297-380G	Human	GST	Sf9 Cells	Full Length
CBL Protein	C272-381G	Human	GST	E.coli	1-375
FANCL Protein	F289-30G	Human	GST	Sf9 Cells	Full Length
Parkin Protein	P215-30G	Human	GST	Sf9 Cells	Full Length
PRPF19 Protein	P270-30G	Human	GST	Sf9 Cells	2-end
STUB1 (CHIP) Protein	C268-30H	Human	HIS	E.coli	Full Length
TRIM32 Protein	T291-30G	Human	GST	Sf9 Cells	2-end

# Ubiquitin Enzymes - E3

Product Name	Cat# No	Species	Tag	Expression System	Sequence
<b>WWP1 Protein</b>	W296-30G	Human	GST	Sf9 Cells	Full Length
<b>VHL/ELOB/ELOC/CUL2/RBX1 complex, Active</b>	U272-380H	Human	HIS	Sf9 Cells	Full length
<b>VHL/ELOB/ELOC/CUL2/RBX1 complex, Active</b>	U272-380H	Human	HIS	Sf9 Cells	Full length

# Ubiquitin Enzymes - DUBs

Product Name	Cat# No	Species	Tag	Expression System	Sequence
AMSH, Active	A565-380H	Human	HIS	Sf9 Cells	Full Length
BAP1, Active	B504-30H	Human	HIS	Sf9 Cells	Full Length
CYLD, Active	C578-380G	Human	GST	Sf9 Cells	1-956
JosD2, Active	I560-380H	Human	HIS	E.coli	Full Length
MINDY1, Protein	M531-30G	Human	GST	Sf9 Cells	Full Length
MINDY4 Protein	M534-30G	Human	GST	Sf9 Cells	Full length
MINDY4, Active	M534-10G	Human	GST	Sf9 Cells	Full length
OTUB1, Protein	O568-30H	Human	HIS	E.coli	Full Length
OTUB2, Active	O569-380H	Human	HIS	E.coli	Full Length
OTUD6B, Protein	O576-30H	Human	HIS	E.coli	Full Length
OTUD7B, Active	O577-381G	Human	GST	Sf9 Cells	1-482
OTULIN, Protein	O585-30H	Human	HIS	E.coli	Full Length
PSMD14 Protein	P501-30G	Human	GST	Sf9 Cells	Full length
SEN1, Active	S531-381H	Human	HIS	E.coli	409-end
SEN2, Active	S532-381H	Human	HIS	E.coli	360-end
SEN6, Active	S536-380H	Human	HIS	E.coli	637-1112
SEN8, Active	S538-380H	Human	HIS	E.coli	Full Length
TNFAIP3 (A20p37), Active	T563-381H	Human	HIS	Sf9 Cells	440-790 a.a.
UCHL1, Active	U522-380H	Human	HIS	E.coli	Full Length
UCHL3, Active	U524-380H	Human	HIS	E.coli	Full Length
UCHL5, Active	U526-380H	Human	HIS	E.coli	Full Length
USP1, Active	U501-31G	Human	GST	Sf9 Cells	Full length
USP10, Active	U510-381G	Human	GST	Sf9 Cells	154-798
USP11, Active	U511-380G	Human	GST	Sf9 Cells	Full Length
USP12, Active	U512-31G	Human	GST	Sf9 Cells	full-length
USP14, Protein	U514-30G	Human	GST	E.coli	Full Length
USP15, Active	U515-380G	Human	GST	Sf9 Cells	Full Length
USP16, Active	U516-381G	Human	GST	Sf9 Cells	187-end
USP2, Active	U502-381H	Human	HIS	E.coli	259-end
USP20, Active	U520-381G	Human	GST	Sf9 Cells	120-end
USP25, Protein	U525-31G	Human	GST	E.coli	162-677 a.a.
USP28, Active	U528-381G	Human	GST	E.coli	155-675a.a.
USP35, Active	U535-381G	Human	GST	Sf9 Cells	395-end
USP46, Protein	U546-30H	Human	HIS	E.coli	Full Length
USP5, Active	U505-380H	Human	HIS	E.coli	Full Length
USP51, Protein	U551-31G	Human	GST	Sf9 Cells	227-end
USP7 (C-Term) Protein	U570-31G	Human	GST	E.coli	560-end

# Ubiquitin Enzymes - DUBs

Product Name	Cat# No	Species	Tag	Expression System	Sequence
USP7 (N-term), Active	U507-381H	Human	HIS	E.coli	25-580a.a.
USP8, Active	U508-381H	Human	HIS	E.coli	630-end
YOD1, Protein	O572-30H	Human	HIS	E.coli	Full Length

## Ubiquitin-like Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
SUMO1 (UBL1) Protein	S293-31H	Human	HIS	E.coli	2-97
SUMO2 Protein	S294-31H	Human	HIS	E.coli	1-93
SUMO3 Protein	S295-31H	Human	HIS	E.coli	1-92
Ubiquitin Protein	U06-54N	Human	TAG-Free	E.coli	1-76
ISG15, Protein	I245-31H	Human	HIS	E.coli	2-157 a. a.
NEDD8, Protein	N263-31H	Human	HIS	E.coli	1-76 a. a.

# Arginine Deiminase & Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
PAD Cocktail, Active	P312-37C		GST	Sf9 Cells	Full Length
PAD1, Active	P312-310G	Human	GST	Sf9 Cells	Full length
PAD2, Active	P312-310BG	Mouse	GST	Sf9 Cells	Full length
PAD2, Active	P312-318BG	Human	GST	Sf9 Cells	Full length
PAD3, Active	P312-310CG	Human	GST	Sf9 Cells	Full length
PAD4, Active	P312-310DG	Human	GST	Sf9 Cells	Full length
PAD6 Protein	P312-30FG	Human	GST	Sf9 Cells	Full Length

## G Proteins

## GTPases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
CDC42, Active	C108-310GH	Human	GST/HIS	E.coli	Full Length
GNAI1, Active	G121-310G	Human	GST	E.coli	2-end
GNAI2, Active	G122-310G	Human	GST	E.coli	2-end
GNAI3, Active	G123-310G	Human	GST	E.coli	2-end
GNAS, Active	G127-310G	Human	GST	Sf9 Cells	Full Length
HRAS (Mature form), Active	R105-310H	Human	HIS	E.coli	2-186
HRAS, Active (Biotinylated)	R105-310HB	Human	Avi-HIS	E.coli	2-186
KRAS (Isoform B), Active	R106B-311H	Human	HIS	E.coli	1-169
KRAS (Mature form), Active	R106-310H	Human	HIS	E.coli	2-186

# GTPases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
KRAS (Mature form, Isoform B), Active	R106B-31H	Human	HIS	E.coli	1-186
KRAS, Active (Biotinylated)	R106-310HB	Human	Avi-HIS	E.coli	2-186
NRAS (Mature form), Active	R107-310H	Human	HIS	E.coli	1-186
NRAS, Active	R107-311H	Human	HIS	E.coli	1-169
NRAS, Active	R107-310AH	Human	Avi-HIS	E.coli	1-186
NRAS, Active (Biotinylated)	R107-31HB	Human	Avi-HIS	E.coli	1-186
NRAS, Active (Biotinylated)	R107-311HB	Human	Avi-HIS	E.coli	1-169
RAC1, Active	R103-310G	Human	GST	E.coli	Full Length
RAC2, Active	R104-310H	Human	HIS	E.coli	Full Length
RAC3, Active	R102-310H	Human	HIS	E.coli	Full Length
RhoA (Mature form), Active	R144-310H	Human	HIS	E.coli	1-190
RhoC (Mature form), Active	R144-310CH	Human	HIS	E.coli	1-190
RhoD (Mature form), Active	R144-310DH	Human	HIS	E.coli	1-207
VCP, Active	V116-310G	Human	GST	Sf9 Cells	Full Length

## Other G Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
GNA11 Protein	G116-30G	Human	GST	E.coli	Full Length
GNA13 Protein	G118-30G	Human	GST	E.coli	Full Length
GNA14 Protein	G119-30G	Human	GST	E.coli	Full Length
GNAZ Protein	G131-30G	Human	GST	E.coli	2-end
HRAS (G12A) Protein	R05-32ATH	Human	HIS	E.coli	2-186
HRAS (G12C) Protein	R05-32CTH	Human	HIS	E.coli	2-186
HRAS (G12S) Protein	R05-32STH	Human	HIS	E.coli	2-186
HRAS (G12V) Protein	R05-32H	Human	HIS	E.coli	2-186
HRAS (G12V) Protein	R05-32VTH	Human	HIS	E.coli	2-186
KRAS (G12C) Protein	R06-32DH	Human	HIS	E.coli	2-186
KRAS (G12C) Protein	R06-32DU	Human	TAG-Free	E.coli	2-186
KRAS (G12C) Protein	R06-32DTH	Human	HIS	E.coli	2-186
KRAS (G12D) Protein	R06-32BH	Human	HIS	E.coli	2-186
KRAS (G12D, Isoform B) Protein	R06B-32BH	Human	HIS	E.coli	1-186
KRAS (G12D, Isoform B) Protein	R06B-321BH	Human	HIS	E.coli	1-169
KRAS (G12V) Protein	R06-32CH	Human	HIS	E.coli	2-186
NRAS (G12C) Protein	R07-32CH	Human	HIS	E.coli	1-169
NRAS (Q61K) Protein	R07-32H	Human	HIS	E.coli	1-186
NRAS (Q61K) Protein	R07-321H	Human	HIS	E.coli	1-169
NRAS (Q61K), Biotinylated	R07-32HB	Human	Avi-HIS	E.coli	1-186
NRAS (Q61K), Biotinylated	R07-321HB	Human	Avi-HIS	E.coli	1-169
RAS Protein	R05-30G	Human	GST	E.coli	Full Length
RGS1 Protein	R39-30H	Human	HIS	E.coli	Full Length
RHEB Protein	R42-30H	Human	HIS	Sf9 Cells	Full Length
RhoA Mutant (S188A) Protein	R44-32AH	Human	HIS	E.coli	1-190

# GAPs/GEFs

Product Name	Cat# No	Species	Tag	Expression System	Sequence
NF1, Active	N120-31H	Human	His	E. coli	1197-1528
RhoGAP, Active	R145-31H	Human	His	E. coli	Full-length
SOS1, Active	S130-31H	Human	His	E. coli	564-1049

# Proteases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
CPB (carboxypeptidase B), Active	C560-31N	Rat	TAG-Free	Yeast cells	109-end
EK (enterokinase), Active	E535-31H	Swine	HIS	Yeast cells	800-end
Immobilized Lys-c, Agarose	L585-31AN	Bacteria	Tag-free	E.coli	206-473 aa
Immobilized Lys-c, Silica	L585-31SN	Bacteria	TAG-Free	E.coli	206-473 aa
Immobilized Thrombin, Agarose	T565-31AN	Human	Native		
Immobilized Trypsin, Agarose	T575-31AN	Swine	TAG-Free	Yeast cells	10-end
Immobilized Trypsin, Silica	T575-31SN	Swine	TAG-Free	Yeast cells	10-end
KEX2 Protease, Active	K525-31N	Yeast	TAG-Free	Yeast cells	110-660 aa
Lys-c (lysyl-endopeptidase), Active	L585-31N	Bacteria	TAG-Free	E.coli	206-473 aa
TEV (Tobacco Etch Virus) Protease, Active	T525-31H	TE Virus	HIS	E.coli	2038-2279 aa
Thrombin, Active	T565-31N	Human	TAG-Free	E.coli	364-end
Trypsin, Active	T575-31N	Swine	TAG-Free	Yeast cells	10-end

# Phosphodiesterases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
PDE10A, Active	P98-31G	Human	GST	Sf9 Cells	71-end
PDE12, Active	P100-310G	Human	GST	Sf9 Cells	Full length
PDE1A, Active	P89-30H	Human	HIS	Sf9 Cells	Full Length
PDE1B, Active	P89-30BG	Human	GST	Sf9 Cells	Full Length
PDE1C, Active	P89-30CG	Human	GST	Sf9 Cells	Full Length
PDE2A, Active	P90-30G	Human	GST	Sf9 Cells	Full Length
PDE3A, Active	P91-31G	Human	GST	Sf9 Cells	669-end
PDE3B, Active	P91-31BG	Human	GST	Sf9 Cells	654-end
PDE4A, Active	P92-31G	Human	GST	Sf9 Cells	332-end
PDE4B, Active	P92-31BG	Human	GST	Sf9 Cells	305-end
PDE4C, Active	P92-31CG	Human	GST	Sf9 Cells	287-end
PDE4D, Active	P92-31DG	Human	GST	Sf9 Cells	225-end
PDE5A, Active	P93-31G	Human	GST	Sf9 Cells	537-end
PDE6A, Active	P94-31G	Human	GST	Sf9 Cells	31-end
PDE6B Protein	P94-34BG	Human	GST	Sf9 Cells	Full length
PDE6C Protein	P94-34CG	Human	GST	Sf9 Cells	Full Length
PDE6D Protein	P94-30DG	Human	GST	Sf9 Cells	Full length
PDE7A, Active	P95-31G	Human	GST	Sf9 Cells	104-end
PDE7B, Active	P95-30BG	Human	GST	Sf9 Cells	Full Length
PDE8A, Active	P96-30G	Human	GST	Sf9 Cells	Full Length
PDE8B, Active	P96-31BG	Human	GST	Sf9 Cells	113-end
PDE9A, Active	P97-30G	Human	GST	Sf9 Cells	Full Length

# Other Enzymes

## Oxidoreductases

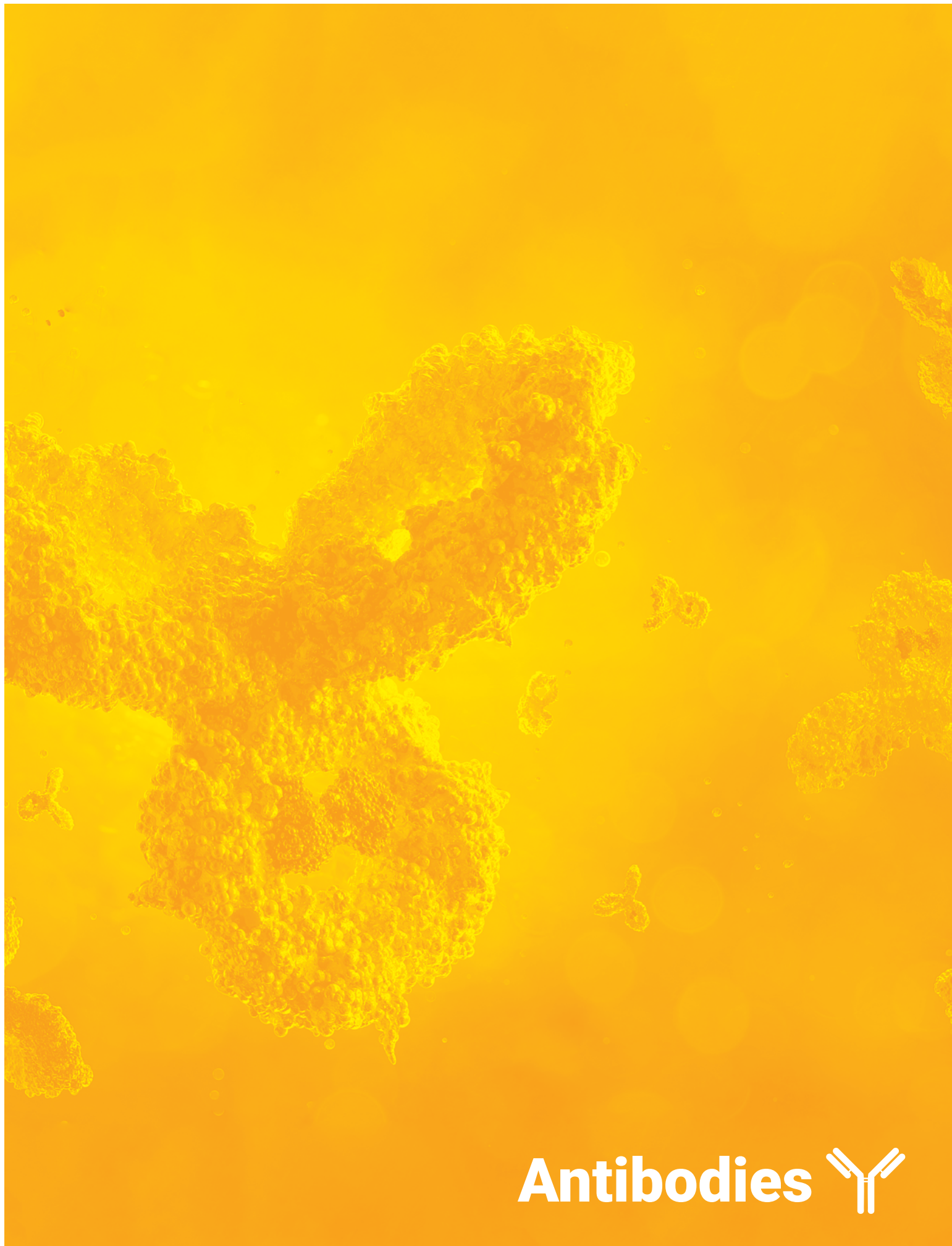
Product Name	Cat# No	Species	Tag	Expression System	Sequence
ALOX12 Protein	A12-34G	Human	GST	Sf9 Cells	Full Length
IDO1 Protein	I211-30G	Human	GST	Sf9 Cells	Full Length
IDO2 Protein	I212-30G	Mouse	GST	Sf9 Cells	Full Length
LOX Protein	L260-31G	Human	GST	Sf9 Cells	169-end
LOXL1 Protein	L261-31G	Human	GST	Sf9 Cells	292-end
LOXL2 Protein	L262-31G	Human	GST	Sf9 Cells	519-end
LOXL3 Protein	L263-31G	Human	GST	Sf9 Cells	526-end
LOXL4 Protein	L264-31G	Human	GST	Sf9 Cells	492-end
SOD1 Protein	S26-30G	Human	GST	E.coli	Full Length
SOD2 Protein	S27-30G	Human	GST	E.coli	Full Length
SOD3 Protein	S28-31G	Human	GST	Sf9 Cells	Full Length
TD02 Protein	T213-30G	Human	GST	Sf9 Cells	Full Length
TET1 Protein	T636-31G	Human	GST	Sf9 Cells	1598-end

## Hydrolases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
CES2, Active	C302-30G	Human	GST	Sf9 Cells	Full length
PLCB1 Protein	P441-30G	Human	GST	Sf9 Cells	Full Length
PLCD4 Protein	P442-30DG	Human	GST	Sf9 Cells	Full Length

## Transferases

Product Name	Cat# No	Species	Tag	Expression System	Sequence
DNA Polymerase $\theta$ , Active	D681-31G	Human	GST	Sf9 Cells	(1-987)
OAS1, Active	O711-30H	Human	HIS	Sf9 Cells	Full length



**Antibodies** 

# Antibodies

## Isoform Specific Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Anti-AKT3	A18-63R	Rabbit	WB	Protein Code: Q9Y243
Anti-AMPK alpha 1	P47-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 483-487 of human AMPK $\alpha$
Anti-CAMK1	C07-60M	Mouse, IgG1, 5H8	WB	Human recombinant full length CAMK1 protein
Anti-CHK2	C48-60M	Mouse, IgG1, 7F9	WB	Human full length recombinant CHK1
Anti-ERK1	M29-60M	Mouse, IgG1, E32	WB, IP	Recombinant full length ERK1 protein
Anti-ERK1	M29-60BM	Mouse, IgG1, E19	WB, IP	Recombinant full length ERK1 protein
Anti-GSK3 alpha	G08-63R	Rabbit	WB	Protein Code P49840
Anti-GSK3 beta	G09-60M	Mouse, IgG1, 6C2	WB	Recombinant full length active GSK3 $\beta$ protein
Anti-HSP90 beta	H36-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 252-256 of human HSP90 $\beta$
Anti-JAK2	J02-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 1005-1009 of human JAK2
Anti-JNK1	M33-63R	Rabbit	WB	Protein Code: P45983
Anti-JNK2	M34-60M	Mouse, IgG2, 2G12	WB, IP	Human recombinant full length JNK2 protein
Anti-MEK1	M02-60M	Mouse, IgG1, 7F1	WB, IP	Recombinant full length humanMEK1 protein
Anti-MEK2	M03-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 392-396 of human MEK2
Anti-p38 alpha	M39-63R	Rabbit	WB	C-terminal peptide of p38 $\alpha$ conjugated to ovalbumin
Anti-p38 beta	M36-63R	Rabbit	WB	C-terminal peptide of p38 $\beta$ conjugated to ovalbumin
Anti-p38 gamma	M37-63R	Rabbit	WB	C-terminal peptide of p38 $\gamma$ conjugated to ovalbumin
Anti-PAK1	P02-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 210-214 of human PAK1
Anti-PKC delta	P64-63CR	Rabbit	WB	Synthesized peptide derived from C-terminus of human PKC $\delta$ protein
Anti-PKC delta	P64-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 643-647 of human PKC $\delta$
Anti-PKC mu	P72-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 736-740 of human PKC $\mu$
Anti-PKC theta	P74-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 674-678 of human PKC $\theta$
Anti-SMAD2	S11-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 218-222 of human SMAD2
Anti-STAT1	S52-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 699-703 of human STAT1
Anti-STAT3	S54-63BR	Rabbit	WB, IP	Synthetic peptide corresponding to amino acids 725-729 of human STAT3
Anti-STAT3	S54-63R	Rabbit	WB	Protein Code: P40763
Anti-STAT4	S55-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 691-695 of human STAT4
Anti-STAT5	S56-63R	Rabbit	WB	Protein Code: P42229

# Isoform Specific Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Anti-STAT5	S56-63BR	Rabbit	WB	Synthetic peptide corresponding to amino acids 692-696 of human STAT5 (protein accession number P42229)

## Tag Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Anti-DYKDDDDK	F51-61M	Mouse, IgG2b, FG4R	WB, ELISA, IP, IS	DYKDDDDK peptide conjugated to KLH
Anti-DYKDDDDK	F51-63R	Rabbit	WB	DYKDDDDK peptide conjugated to KLH
Anti-GFP	G46-60M	Mouse, IgG1, GF28R	WB	GFP N-terminal peptide with KLH conjugate from Aequorea Victoria Jellyfish
Anti-GST	G52-61M	Mouse, IgG2a, G2R	WB	Purified GST
Anti-HA	H98-61M	Mouse, IgG3, C10	WB, ELISA, IP, IS	
Anti-HA	H98-63R	Rabbit Polyclonal	WB, ELISA, IP, IS	
Anti-His	H99-61M	Mouse, IgG2b, EH158	WB, ELISA, IP, IS	HHHHHH peptide conjugated to KLH
Anti-His	H99-63R	Rabbit	WB, ELISA, IP, IS	
Anti-Myc	M86-61M	Mouse, IgG1, M19	WB, ELISA, IP, IS	EQKLISEED peptide conjugated to KLH
Anti-RFP	R46-61M	Mouse, IgG1, RF5R	WB	RFP from the Discosoma sea anemone N-terminal peptide-KLH conjugates
Anti-V5	V05-61M	Mouse, IgG1, V4RR	WB, ELISA, IP, IS	GKPIPPLGLDST peptide conjugated to KLH

## Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Alpha Synuclein Antibody	S22-62R	Rabbit	WB, IHC, ELISA	Recombinant Human alpha Synuclein Fibril
Alpha Synuclein Antibody (pSer129)	S22-635R	Rabbit	WB, ICC/IF, IHC, ELISA	Synthetic peptide of Human Alpha Synuclein pSer129 (40-140aa), conjugated to Keyhole Limpet Haemocyanin (KLH).
Alpha Synuclein Antibody, 10H7	S22-61BM	Mouse / IgG1 / 10H7	WB, DB, ICC/IF, ELISA	Mouse alpha synuclein aggregate
Alpha Synuclein Antibody, 3C11	S22-61M	Mouse / IgG1 / 3C11	WB, DB, ICC/IF, ELISA	Human alpha synuclein monomer
Alpha Synuclein Antibody, 3F8	S22-61CM	Mouse / IgG1 / 3F8	WB, DB, 1HC, ICC/IF, ELISA	Mouse alpha synuclein aggregate
Alpha Synuclein Antibody, 4F1	S22-61DM	Mouse / IgG1 / 4F1	WB, DB, IHC, ICC/IF, ELISA	Mouse alpha synuclein aggregate
Anti-4EBP1	E35-563R	Rabbit, IgG	WB, ELISA, ICC	4EBP1 antibody was raised against a 14 amino acid synthetic peptide from near the carboxy terminus of human 4EBP1
Anti-ADD1	A38-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 724-728 of human ADD1 (P35611)
Anti-AFX	A42-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 195-199 of human AFX
Anti-AKT (pan)	A16-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 471-475 of human AKT

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-AKT (pan)</b>	A16-63BR	Rabbit	WB	Synthetic peptide corresponding to amino acids 313-317 of human AKT1 (protein accession number P31749)
<b>Anti-AKT1</b>	A16-60M	Mouse, Monoclonal, 9A4	WB	Recombinant full length active AKT1 protein
<b>Anti-AKT1</b>	A16-63CR	Rabbit, IgG	WB, ELISA, IF, IHC	AKT1 antibody was raised against a 16 amino acid synthetic peptide from near the amino-terminus of human AKT1
<b>Anti-AKT2</b>	A17-61A	Rat, IgG2a	ELISA, IHC, WB	The antibody was produced against synthesized peptide corresponding to internal residues of the human AKT2 protein.
<b>Anti-AKT2/3</b>	A17-63BR	Rabbit	WB	Protein Code: P31751
<b>Anti-AKT3</b>	A18-61M	Mouse, IgG1	ELISA, IHC, WB	The antibody was produced against synthesized peptide corresponding to internal residues of the human AKT3 protein.
<b>Anti-ALK2</b>	A06-63R	Rabbit, IgG	WB, ELISA	ALK2 antibody was raised against a 14 amino acid synthetic peptide near the amino terminus of the human ALK2
<b>Anti-ALK4</b>	A07-63R	Rabbit, IgG	WB, ELISA	ALK4 antibody was raised against a 14 amino acid synthetic peptide near the carboxy terminus of the human ALK4
<b>Anti-ALK7</b>	A08-63R	Rabbit, IgG	WB, ELISA	ALK7 antibody was raised against a 15 amino acid synthetic peptide near the amino terminus of the human ALK7
<b>Anti-alpha 1-Actin</b>	A03-360M	Mouse, IgG1, C4	WB, IF, IHC	Chicken gizzard actin
<b>Anti-AMPK alpha 1/2</b>	P47-63BR	Rabbit	WB	Synthetic peptide corresponding to amino acids 170-174 of human AMPK
<b>Anti-Amyloid beta</b>	A06-563R	Rabbit, IgG	WB, ELISA, IF, IHC	Amyloid beta antibody was raised against a synthetic peptide corresponding to the amino terminus (aa 1-14) of human beta amyloid conjugated to KLH using maleimide.
<b>Anti-AOS1</b>	A80-63R	Rabbit	WB	Synthetic peptide of c-terminal amino acid sequence of human AOS1 (Q9UBE0)
<b>Anti-APBA1</b>	A86-363R	Rabbit, IgG	WB	Synthetic peptide corresponding to amino acid residues from the C-terminal region conjugated to KLH
<b>Anti-APC1</b>	A87-363R	Rabbit, IgG	WB, ELISA, IHC	APC1 antibody was raised against a 17 amino acid synthetic peptide near the carboxy terminus of human APC1
<b>Anti-APC1 (Ser355)</b>	A87-363BR	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to amino acids 351-359 of Human APC1.
<b>Anti-APC10</b>	A95-363R	Rabbit, IgG	WB, ELISA, IF, IHC	APC10 antibody was raised against a 16 amino acid synthetic peptide near the center of human APC10
<b>Anti-APC11</b>	A96-363R	Rabbit, IgG	WB, ELISA, IF	APC11 antibody was raised against an 18 amino acid synthetic peptide near the center of human APC11
<b>Anti-APC13</b>	A97-363R	Rabbit, IgG	WB, ELISA, IF	APC13 antibody was raised against a 17 amino acid synthetic peptide near the center of human APC13
<b>Anti-APC2</b>	A88-363R	Rabbit, IgG	WB, ELISA, ICC, IF	APC2 antibody was raised against an 18 amino acid synthetic peptide near the center of human APC2

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-APC3</b>	A89-363R	Rabbit, IgG	WB, ELISA, IF, IHC	APC3 antibody was raised against a 19 amino acid synthetic peptide near the carboxy terminus of human APC3
<b>Anti-APC3 (CDC27)</b>	A89-363BR	Rabbit, IgG	ELISA, IF, WB	The antibody was produced against synthesized peptide corresponding to amino acids 422-430 of human Cdc27
<b>Anti-APC4</b>	A90-363R	Rabbit, IgG	WB, ELISA, IHC	APC4 antibody was raised against a 14 amino acid synthetic peptide near the amino terminus of human APC4
<b>Anti-APC5</b>	A91-363R	Rabbit, IgG	WB, ELISA, IF, IHC	APC5 antibody was raised against a 17 amino acid synthetic peptide near the center of human APC5
<b>Anti-APC6</b>	A92-363R	Rabbit, IgG	WB, ELISA, IF, IHC	APC6 antibody was raised against an 18 amino acid synthetic peptide near the amino terminus of human APC6
<b>Anti-APC7</b>	A93-363R	Rabbit, IgG	WB, ELISA, IF	APC7 antibody was raised against an 18 amino acid synthetic peptide near the carboxy terminus of human APC7
<b>Anti-APC8</b>	A94-363R	Rabbit, IgG	WB, ELISA, ICC	APC8 antibody was raised against a 17 amino acid synthetic peptide near the center of human APC8
<b>Anti-ASK1</b>	M13-60M	Mouse, A1	WB	Human recombinant 649-946 ASK1
<b>Anti-ASK1</b>	M13-60BM	Mouse, A23	WB	Human recombinant 649-946 ASK1
<b>Anti-ASK1</b>	M13-63BR	Rabbit, IgG	WB, ELISA, ICC, IF	ASK1 antibody was raised against a 16 amino acid synthetic peptide from near the carboxy terminus of human ASK1.
<b>Anti-ASK1</b>	M13-63R	Rabbit, IgG	WB, ELISA, ICC, IF	ASK1 antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human ASK1. This sequence is different from that of mouse by last two amino acids.
<b>Anti-ASK1</b>	M13-63CR	Rabbit, IgG	WB, ELISA	ASK1 antibody was raised against a 16 amino acid synthetic peptide from near the amino terminus of human ASK1
<b>Anti-ATF2</b>	A10-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 60-64 of human ATF2
<b>Anti-ATG13</b>	A113-363R	Rabbit, IgG	WB, ELISA, IF, IHC	ATG13 antibody was raised against a 15 amino acid synthetic peptide near the carboxy terminus of human ATG13
<b>Anti-ATG5</b>	A105-363R	Rabbit, IgG	WB, ELISA, IF, IHC	ATG5 antibody was raised against a 16 amino acid synthetic peptide from near the amino terminus of human ATG5.
<b>Anti-AURORA B</b>	A31-63R	Rabbit, IgG	ELISA, IF, WB	The antibody was produced against synthesized peptide corresponding to amino acids 277-238 of Human Aurora Kinase B protein.
<b>Anti-AXL</b>	A34-63R	Rabbit, IgG	WB, ELISA	Peptide at the N-terminal residues of human AXL protein.
<b>Anti-BAD</b>	B03-363R	Rabbit, IgG	WB, ELISA, IF, IHC	BAD antibody was raised against a peptide corresponding to 15 amino acids near the C-terminus of human BAD
<b>Anti-BAG1</b>	B05-363BR	Rabbit, IgG	WB, ELISA, IF, IHC	BAG1 antibody was raised against a 14 amino acid synthetic peptide from near the amino terminus of human BAG1

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-BAG1</b>	B05-363R	Rabbit, IgG	WB, ELISA	BAG1 antibody was raised against a 14 amino acid synthetic peptide from near the carboxy terminus of human BAG1
<b>Anti-BAK1</b>	B09-363R	Rabbit, IgG	WB, ELISA, ICC	BAK1 antibody was raised against a peptide corresponding to 13 amino acids near the amino-terminus of human BAK1
<b>Anti-BAX</b>	B14-363R	Rabbit, IgG	WB, ELISA, ICC, IF	BAX antibody was raised against a peptide corresponding to 16 amino acids near the amino-terminus of human BAX
<b>Anti-BCL2</b>	B25-63R	Rabbit	WB, IP	Synthetic peptide corresponding to amino acids 54-58 of human BCL2
<b>Anti-BCL2</b>	B25-63BR	Rabbit, IgG	WB, ELISA, IF, IHC	BCL2 antibody was raised against a peptide corresponding to 15 amino acids near the N terminus of human BCL2
<b>Anti-BCL2</b>	B25-63CR	Rabbit, IgG	WB, ELISA, ICC	BCL2 antibody was raised against a peptide corresponding to 14 amino acids near the middle of human BCL2
<b>Anti-BCL2L1</b>	B26-363R	Rabbit, IgG	WB, ELISA	BCL2L1 antibody was raised against a peptide corresponding to 14 amino acids near the carboxy terminus of human BCL2L1
<b>Anti-BCL2L10</b>	B26-363JR	Rabbit, IgG	WB, ELISA, ICC	BCL2L10 antibody was raised against a 16 amino acid synthetic peptide from near the amino terminus of human BCL2L10
<b>Anti-BCL2L14</b>	B26-363NR	Rabbit, IgG	WB, ELISA, IF, IHC	BCL2L14 antibody was raised against a peptide corresponding to 15 amino acids near the C-terminus of human BCL2L14
<b>Anti-beta-Actin</b>	A03-63BR	Rabbit	WB	Protein Code: P60709
<b>Anti-beta-Actin</b>	A03-60BM	Mouse, IgG2b, BA3R	WB	Slightly modified human $\beta$ -cytoplasmic actin (2-15aa) peptide
<b>Anti-BID</b>	B59-363BR	Rabbit, IgG	WB, ELISA	BID antibody was raised against a peptide corresponding to 14 amino acids near the carboxy terminus of human BID
<b>Anti-BID</b>	B59-363R	Rabbit, IgG	WB, ELISA, IF, IHC	BID antibody was raised against a peptide corresponding to 14 amino acids near the carboxy terminus of human BID
<b>Anti-BIRC2</b>	B279-63R	Rabbit, IgG	WB, ELISA, IF, IHC	BIRC2 antibody was raised against a synthetic peptide corresponding to 14 amino acids at the C-terminus of human BIRC2
<b>Anti-BMPR2</b>	B06-63R	Rabbit, IgG	WB, ELISA, IHC	Recombinant human BMPR2 protein corresponding to amino acids 295-552
<b>Anti-BOK</b>	B70-363R	Rabbit, IgG	ELISA, IHC	BOK antibody was raised against a 16 amino acid synthetic peptide near the amino terminus of human BOK
<b>Anti-BRAF</b>	B08-63R	Rabbit, IgG	WB, ELISA	BRAF antibody was raised against an 18 amino acid synthetic peptide near the center of human BRAF
<b>Anti-BRSK1</b>	B13-63BR	Rabbit, IgG	WB, ELISA, IF, IHC	BRSK1 antibody was raised against a 28 amino acid synthetic peptide from near the center of human BRSK1
<b>Anti-BRSK1</b>	B13-63R	Rabbit, IgG	WB, ELISA, IF, IHC	BRSK1 antibody was raised against a 28 amino acid synthetic peptide from near the carboxy terminus of human BRSK1
<b>Anti-BRSK2</b>	B14-63R	Rabbit, IgG	WB, ELISA, IF, IHC	BRSK2 antibody was raised against a 23 amino acid synthetic peptide from near the carboxy terminus of human BRSK2.

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-BTK</b>	B10-63R	Rabbit, IgG	WB, ELISA, ICC, IF	BTK antibody was raised against a synthetic peptide corresponding to 16 amino acids near the amino-terminus of human BTK
<b>Anti-BUB1</b>	B11-63BR	Rabbit, IgG	WB, ELISA, ICC, IF	BUB1 antibody was raised against a 22 amino acid synthetic peptide from near the carboxy terminus of human BUB1
<b>Anti-BUB1</b>	B11-63R	Rabbit, IgG	WB, ELISA, ICC, IF	BUB1 antibody was raised against a 19 amino acid synthetic peptide from near the amino terminus of human BUB1
<b>Anti-BUB1B</b>	B12-60M	Mouse, IgG1	ELISA, IF, IP, WB	The antibody was produced against synthesized peptide corresponding to amino acid residues 1-350 of human BUBR1 Kinase protein.
<b>Anti-c-JUN</b>	C05-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 71-75 of human c-JUN
<b>Anti-c-KIT</b>	K06-63R	Rabbit, IgG	ELISA, IHC	Rabbit polyclonal c-KIT antibody was raised against a 16 amino acid peptide from near the C-terminal residues of human c-kit protein.
<b>Anti-CAMK2</b>	C11-63R	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to amino acids 6-23 of Human CaM Kinase II protein.
<b>Anti-CAMK4</b>	C15-63R	Rabbit, Antiserum	ELISA, WB	The antibody was produced against synthesized peptide corresponding to amino acids 305-323 of Human CaM Kinase IV protein.
<b>Anti-CASK</b>	C19-63R	Rabbit, IgG	WB, ELISA, IF	CASK antibody was raised against a 16 amino acid synthetic peptide near the amino terminus of human CASK
<b>Anti-CASKIN1</b>	C13-363R	Rabbit, IgG	WB, ELISA	CASKIN1 antibody was raised against a 17 amino acid synthetic peptide near the center of human CASKIN1
<b>Anti-Caspase-3</b>	C603-363CR	Rabbit, IgG	ELISA, IHC	Caspase-3 antibody was raised against a 17 amino acid synthetic peptide near the center of human Caspase-3.
<b>Anti-CD5L (CT-2)</b>	C61-363BR	Rabbit, IgG	WB, ELISA, IF, IHC	CD5L (CT-2) antibody was raised against a 16 amino acid synthetic peptide near the amino terminus of human CD5L (CT-2)
<b>Anti-CD5L (CT-2)</b>	C61-363R	Rabbit, IgG	WB, ELISA, IHC	CD5L (CT-2) antibody was raised against a 13 amino acid synthetic peptide near the carboxy terminus of human CD5L (CT-2)
<b>Anti-CDC25A</b>	C04-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 74-78 of human CDC25A
<b>Anti-CDC42</b>	C08-363R	Rabbit, IgG	WB, ELISA	CDC42 antibody was raised against a 17 amino acid synthetic peptide near the amino terminus of human CDC42
<b>Anti-CDK1</b>	C22-63R	Rabbit, Antiserum	ELISA, IHC, IP, WB	The antibody was produced against synthesized peptide corresponding to the C-terminus of the human protein conjugated to KLH.
<b>Anti-CDK2</b>	C29-63R	Rabbit, Antiserum	ELISA, IHC, IP, WB	The antibody was produced against synthesized peptide corresponding to the C-Terminus of the human CDK2 protein conjugated to KLH.

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-CDK4</b>	C31-63R	Rabbit, Antiserum	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to the C-Terminus of the human CDK4 (p34) protein conjugated to KLH.
<b>Anti-CDK5</b>	C33-60M	Mouse, IgG1, 1H3	WB, IHC	Purified rat Cdk5
<b>Anti-CDK5</b>	C33-63R	Rabbit, Antiserum	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to the C-Terminus of human CDK5 protein conjugated to KLH.
<b>Anti-CDK9</b>	C40-63R	Rabbit, Antiserum	ELISA, IHC, IP, WB	The antibody was produced against synthesized peptides corresponding to the C-terminal and N-terminal domains of human CDK9 (PITALRE) protein.
<b>Anti-CHK2</b>	C48-63R	Rabbit, IgG	WB, ELISA, ICC	CHK2 antibody was raised against a synthetic peptide corresponding to amino acids near the amino terminus of human CHK2
<b>Anti-Choline Acetyltransferase</b>	C309-363G	Goat, IgG	WB, IHC	Native choline acetyltransferase purified from human placenta
<b>Anti-CREB</b>	C49-63R	Rabbit	WB, IP	Synthetic peptide corresponding to amino acids 699-703 of human STAT1
<b>Anti-Cyclin A</b>	C78-363R	Rabbit, IgG	WB, ELISA, IP	Anti-Cyclin A Antibody was produced by repeated immunizations with a recombinant protein corresponding to the human cyclin A.
<b>Anti-Cyclin A</b>	C78-363BR	Rabbit, Antiserum	ELISA, IP, WB	The antibody was produced against a recombinant protein corresponding to the human Cyclin A protein.
<b>Anti-Cyclin B1</b>	C79-362R	Rabbit, Antiserum	ELISA, IP, WB	The antibody was produced against full length fusion protein corresponding to the human Cyclin B1 protein.
<b>Anti-Cyclin D1</b>	C81-362R	Rabbit, Antiserum	ELISA, IP, WB	The antibody was produced against full length fusion protein corresponding to the human Cyclin D1 gene sequence.
<b>Anti-Cyclin E</b>	C82-363R	Rabbit, Antiserum	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to a region near the carboxy terminus of human cyclin E conjugated to KLH.
<b>Anti-Cyclin E2</b>	C82-363BR	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to amino acids at the carboxyl terminus of Cyclin E2 protein
<b>Anti-Cyclin O</b>	C85-363BR	Rabbit, IgG	WB, ELISA	CyclinO antibody was raised against a 13 amino acid synthetic peptide from near the center of human CyclinO
<b>Anti-Cyclin O</b>	C85-363R	Rabbit, IgG	WB, ELISA	CyclinO antibody was raised against a 14 amino acid synthetic peptide from near the amino terminus of human CyclinO.
<b>Anti-CyclinT1</b>	C86-363BR	Rabbit, IgG	WB, ELISA	CyclinT1 antibody was raised against an 18 amino acid peptide near the carboxy terminus of human CyclinT1
<b>Anti-Cytokeratin 5 (KRT5)</b>	C91-563R	Rabbit, IgG	WB, ELISA	Synthetic peptide conjugated to KLH

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Anti-DAPK2	D02-63R	Rabbit, IgG	WB, ELISA, IHC	DAPK2 antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human DAPK2. This antigenic peptide is identical to the corresponding amino acids of mouse origin.
Anti-DAPK3	D03-63R	Rabbit, IgG	WB, ELISA, ICC, IF	DAPK3 antibody was raised against a peptide corresponding to amino acids near the center of human DAPK3.
Anti-DCAMKL1	D14-63R	Rabbit, IgG	WB, ELISA, IF, IHC	DCAMKL1 antibody was raised against a 14 amino acid synthetic peptide near the amino terminus of human DCAMKL1
Anti-DCAMKL2	D15-63R	Rabbit, IgG	WB, ELISA, IHC	DCAMKL2 antibody was raised against a 14 amino acid synthetic peptide near the carboxy terminus of human DCAMKL2.
Anti-DCAMKL3	D16-63R	Rabbit, IgG	WB, ELISA, IF	DCAMKL3 antibody was raised against a 14 amino acid synthetic peptide near the amino terminus of human DCAMKL3
Anti-DLK1	M20-63R	Rabbit, IgG	WB, ELISA	DLK1 antibody was raised against an 18 amino acid synthetic peptide near the carboxy terminus of human DLK1
Anti-DNMT3L	D353-63R	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to amino acids 152-164 of human DNMT3L (DNA (cytosine-5-)-methyltransferase-3-like).
Anti-DOK1	D42-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 360-364 of human DOK1 (Q99704)
Anti-DRAK2 (STK17B)	S34-63R	Rabbit, IgG	WB, ELISA, ICC	DRAK2 (STK17B) antibody was raised against a peptide corresponding to amino acids 351 to 365 of human DRAK2 (STK17B)
Anti-DYRK1A	D09-63R	Rabbit, IgG	WB, ELISA, IF	DYRK1A antibody was raised against a 17 amino acid synthetic peptide near the carboxy terminus of human DYRK1A.
Anti-DYRK2	D10-63R	Rabbit, IgG	WB, ELISA, ICC, IF	DYRK2 antibody was raised against an 18 amino acid synthetic peptide near the carboxy terminus of human DYRK2
Anti-EGFR	E10-63BR	Rabbit	WB	Synthetic peptide corresponding to amino acids 1195- 1199 of human EGFR (protein accession number P00533)
Anti-EGFR	E10-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 1090- 1094 of human EGFR (protein accession number P00533)
Anti-EIF2AK2 (PKR)	P80-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 449-453 of human EIF2AK2
Anti-EIF2AK2 (PKR)	P80-63BR	Rabbit, IgG	WB, ELISA, IF, IHC	EIF2AK2 antibody was raised against a peptide corresponding to 14 amino acids near the carboxy terminus of human EIF2AK2
Anti-EIF2AK2 (PKR)	P80-63CR	Rabbit, IgG	WB, ELISA, IHC	EIF2AK2 antibody was raised against a peptide corresponding to 14 amino acids near the amino terminus of human EIF2AK2
Anti-eIF4E	E34-63R	Rabbit	WB	Synthesized peptide corresponds to amino acid 209 of human eIF4E protein.
Anti-ELK1	E64-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 387-391 of human ELK1

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Anti-ERK1	M29-60CM	Mouse, IgG2b, E40	IP	Recombinant full length ERK1 protein
Anti-ERK1/2	M29-60EM	Mouse, IgG1, E31R	WB	Human full length recombinant ERK1
Anti-ERK2	M28-63R	Rabbit	WB	Human ERK2 c-terminal peptide-KLH conjugates
Anti-EZH1	E95-363R	Rabbit, IgG	WB, ELISA, IF, IHC	EZH1 antibody was raised against an 18 amino acid synthetic peptide near the amino terminus of human EZH1
Anti-EZH2	E96-363R	Rabbit, IgG	WB, ELISA, ICC, IF	EZH2 antibody was raised against a 17 amino acid synthetic peptide near the amino terminus of human EZH2.
Anti-FGF4	F824-63R	Rabbit, IgG	WB, ELISA, ICC, IF	FGF4 antibody was raised against a 18 amino acid synthetic peptide near the carboxy terminus of the human FGF4
Anti-FLT3	F12-63R	Rabbit, IgG	WB, ELISA, IF	FLT3 antibody was raised against an 18 amino acid synthetic peptide near the center of human FLT3
Anti-FOXO1	F19-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 317-321of human FOXO1
Anti-FOXO3	F70-363R	Rabbit, IgG	WB, ELISA, ICC, IF	Anti-FOXO3 antibody was raised against an 18 amino acid synthetic peptide near the amino terminus of human FOXO3.
Anti-GAB1	G03-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 625-629 of human GAB1
Anti-GAPDH	G13-61M	Mouse, IgG1, GA1R	WB, ELISA	GAPDH from rabbit
Anti-GAPDH	G13-63R	Rabbit	WB	Synthetic peptide corresponding to a portion of the human G3PDH
Anti-GATA1	G19-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 308-312 of human GATA1
Anti-GFAP	G28-563R	Rabbit, IgG	WB, IF	Recombinant and purified bovine GFAP.
Anti-GLUT1	G837-363R	Rabbit, IgG	WB, ELISA	Synthetic peptide conjugated to KLH
Anti-GSK3 alpha/beta	G09-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 214-218 of human GSK3β
Anti-HDAC1	H83-63R	Rabbit	WB	Protein Code Q13547
Anti-HDAC11	H93-63R	Rabbit	WB	Synthetic peptide corresponding to c-terminus of human HDAC11 conjugated to KLH.
Anti-HDAC2	H84-63R	Rabbit	WB	Protein Code Q92769
Anti-HDAC4	H86-63R	Rabbit	WB	Protein Code P56524
Anti-HDAC4	H86-63BR	Rabbit, IgG	ELISA, IF, IC, WB	The antibody was produced against synthesized peptide corresponding to the internal region surrounding 600-650 aa of human HDC4 protein
Anti-HDAC5	H87-63R	Rabbit	WB	Synthetic peptide corresponding to c-terminus of human HDAC5 conjugated to KLH.
Anti-HDAC5	H87-63BR	Rabbit, IgG	ELISA, IHC, IF, WB	The antibody was produced against synthesized peptide corresponding to the internal region surrounding phosphoserine 661 of human HDAC5
Anti-HDAC6	H88-63R	Rabbit	WB	Synthetic peptide corresponding to c-terminus of human HDAC6 conjugated to KLH.

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Anti-HDAC7	H89-63R	Rabbit	WB	Synthetic peptide corresponding to c-terminus of human HDAC7 conjugated to KLH.
Anti-HDAC8	H90-63R	Rabbit	WB	Protein Code Q9BY41
Anti-HDAC9	H91-63R	Rabbit	WB	Protein Code Q9UKV0
Anti-HER2	E27-63R	Rabbit, IgG	WB, ELISA, IF, IHC	HER2 antibody was raised against a 19 amino acid synthetic peptide near the carboxy terminus of human HER2
Anti-HER3	E28-62R	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against HER3 fusion protein corresponding to amino acids 1283 and 1323 at the C-terminus of human HER3
Anti-Histone H3	H12-63R	Rabbit, antiserum	ELISA, WB	The antibody was produced against synthesized peptide corresponding to the C-Terminus region of human histone 3
Anti-HSF1	H25-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 305-309 of human HSF1
Anti-HSP27	H30-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 80-84 of human HSP27 (protein accession number P04792)
Anti-HSP27	H30-60M	Mouse, IgG1	ELISA, WB	The antibody was produced against prokaryotic recombinant protein corresponding to the full length human HSP27 protein
Anti-IKK alpha	C51-63BR	Rabbit, IgG	WB, ELISA, ICC, IF	IKK alpha antibody was raised against a peptide corresponding to amino acids 699 to 715 of human IKK alpha, which differs from corresponding murine sequence by one amino acid.
Anti-IKK alpha	C51-63CR	Rabbit, IgG	WB, ELISA, ICC	IKK alpha antibody was raised against a peptide corresponding to amino acids 716 to 734 of human IKK alpha, which differs from corresponding murine sequence by four amino acids.
Anti-IKK alpha	C51-63R	Rabbit, IgG	WB, ELISA, ICC, IF	IKK alpha antibody was raised against a peptide corresponding to amino acids 716 to 734 of human IKK alpha, which differs from corresponding murine sequence by four amino acids.
Anti-IKK beta	I03-63R	Rabbit, IgG	WB, ELISA, ICC, IF	IKK beta antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human IKK beta, which differs from corresponding murine sequence by one amino acid.
Anti-IKK epsilon	I04-63R	Rabbit, IgG	WB, ELISA, IHC	IKK epsilon antibody was raised against a peptide corresponding to amino acids 701 to 716 of human IKK epsilon
Anti-IKK gamma	I25-63R	Rabbit, IgG	WB, ELISA, ICC	IKK gamma antibody was raised against a peptide corresponding to amino acids 400 to 416 of human IKK gamma, which are identical to those of mouse homologue.
Anti-IL-21	I881-463R	Rabbit, IgG	WB, ELISA, ICC	Anti-IL-21 antibody was raised against a synthetic peptide corresponding to 15 amino acids near the center of human IL-21 precursor.

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-IRAK1</b>	I09-63R	Rabbit, IgG	WB, ELISA, ICC, IF, IP	IRAK1 antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human IRAK1
<b>Anti-IRAK2</b>	I10-63BR	Rabbit, IgG	WB, ELISA, ICC, IF	IRAK2 antibody was raised against a peptide corresponding to 17 amino acids near the carboxy terminus of mouse IRAK2a which is common to all four isoforms
<b>Anti-IRAK2</b>	I10-63CR	Rabbit, IgG	WB, ELISA	IRAK2 antibody was raised against a peptide corresponding to 16 amino acids near the carboxy terminus of mouse IRAK2a which is common to all four isoforms
<b>Anti-IRAK2</b>	I10-63R	Rabbit, IgG	WB, ELISA, ICC, IF	IRAK2 antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human IRAK2
<b>Anti-IRAK3</b>	I11-63R	Rabbit, IgG	WB, ELISA, IF, IHC	IRAK3 antibody was raised against a peptide corresponding to amino acids from the carboxy terminus of human IRAK3
<b>Anti-IRAK4</b>	I12-63R	Rabbit, IgG	WB, ELISA, ICC, IF	IRAK4 antibody was raised against a synthetic peptide corresponding to amino acids at the carboxy terminus of human IRAK4
<b>Anti-IRS1</b>	I40-363R	Rabbit, IgG	WB, ELISA, ICC, IF	IRS1 antibody was raised against a 16 amino acid synthetic peptide from near the center of human IRS1
<b>Anti-JMJD1C</b>	J48-363R	Rabbit, IgG	WB, ELISA, IF, IHC	JMJD1C antibody was raised against a 20 amino acid synthetic peptide from near the amino terminus of human JMJD1C
<b>Anti-JNK2/3</b>	M34-63R	Rabbit	WB	C-terminal peptide of JNK3 conjugated to ovalbumin
<b>Anti-KDM1A</b>	K421-63BR	Rabbit, IgG	WB, ELISA, IF, IHC	KDM1A antibody was raised against a 16 amino acid synthetic peptide from near the center of human KDM1A
<b>Anti-KDM1A</b>	K421-63R	Rabbit, IgG	WB, ELISA	KDM1A antibody was raised against a 17 amino acid synthetic peptide from near the amino terminus of human KDM1A
<b>Anti-KDM3A</b>	K423-63R	Rabbit, IgG	WB, ELISA, IF, IHC	KDM3A antibody was raised against a 16 amino acid synthetic peptide from near the amino terminus of human KDM3A
<b>Anti-KDM3B</b>	K423-63BR	Rabbit, IgG	WB, ELISA, IF, IHC	KDM3B antibody was raised against a 20 amino acid synthetic peptide from near the center of human KDM3B
<b>Anti-KDM4A</b>	K424-63R	Rabbit, IgG	WB, ELISA, IHC	KDM4A antibody was raised against a 14 amino acid synthetic peptide from near the center of human KDM4A
<b>Anti-KDM4B</b>	K424-63BR	Rabbit, IgG	WB, ELISA	KDM4B antibody was raised against a 16 amino acid synthetic peptide from near the center of human KDM4B
<b>Anti-KDM6B</b>	K426-63BR	Rabbit, IgG	WB, ELISA, ICC, IF	KDM6B antibody was raised against a 16 amino acid synthetic peptide from near the amino terminus of human KDM6B
<b>Anti-KDM8</b>	K428-63R	Rabbit, IgG	WB, ELISA, IF, IHC	KDM8 antibody was raised against an 18 amino acid synthetic peptide from near the carboxy terminus of human KDM8

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-LIMK1</b>	L04-63R	Rabbit, IgG	ELISA, IF, WB	The antibody was produced against synthesized peptide corresponding to amino acids 630-647 of human LIM Kinase protein.
<b>Anti-LKB1 (STK11)</b>	L15-63R	Rabbit, IgG	WB, ELISA	LKB1 antibody was raised against a 15 amino acid synthetic peptide from near the carboxy terminus of human LKB1
<b>Anti-MAP2</b>	M110-563R	Rabbit, IgG	ELISA, IHC	MAP2 antibody was raised against a recombinant human MAP2 protein 339-554aa
<b>Anti-MAPKAPK2/3</b>	M40-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 332-336 of human MAPKAPK2
<b>Anti-MARCKS</b>	M02-562R	Rabbit, IgG	WB, IF	Recombinant protein expressed in and purified from E. coli
<b>Anti-MBP</b>	M42-563C	Chicken, IgY	WB, IF	Three peptide sequences conserved in higher vertebrate MBP proteins
<b>Anti-MCAM</b>	M141-563R	Rabbit, IgG	WB, ELISA	N/A
<b>Anti-MEK1/2</b>	M02-63R	Rabbit	WB	Synthesized peptide corresponding to human MEK1 (210-220 aa) protein
<b>Anti-mGLUR5</b>	G834-463R	Rabbit, IgG	WB, IHC	Peptide from the C-terminal region of rat mGLUR5 and rat mGLUR1a.
<b>Anti-mTOR</b>	F17-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 2446-2450 of human mTOR
<b>Anti-mTOR</b>	F17-63BR	Rabbit, IgG	WB, ELISA, ICC	mTOR antibody was raised against a 15 amino acid synthetic peptide from near the amino terminus of human mTOR
<b>Anti-MYT1L</b>	M92-363R	Rabbit, IgG	WB, ELISA, IF	Rabbit polyclonal MYT1L antibody was raised against an 18 amino acid peptide near the carboxy terminus of human MYT1L
<b>Anti-NEK2</b>	N03-63R	Rabbit, IgG	ELISA, IF, WB	The antibody was produced against synthesized peptide corresponding to amino acids 287-299 of human NEK2 protein
<b>Anti-Nestin (NES)</b>	N122-561M	Mouse, IgG	ELISA, IHC	Recombinant human Nestin expressed in bacteria and highly purified. The specific region was amino acids 317 to 630 of the human protein. This is a region which is 55% identical to rat Nestin and 57% identical to mouse Nestin.
<b>Anti-Neurofilament-L (NF-L)</b>	N125-561M	Mouse, IgG, DA2	WB, IF	Preparation of porcine intermediate filaments
<b>Anti-NF-κB p105</b>	N12-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 930-934 of human NF-κB p105
<b>Anti-NIK</b>	N22-63R	Rabbit, IgG	WB, ELISA	NIK antibody was raised against a peptide corresponding to amino acids 931 to 947 of human NIK
<b>Anti-NOTCH 2</b>	N42-363R	Rabbit, IgG	WB, ELISA, IHC	This whole rabbit serum was prepared by repeated immunizations with a synthetic peptide corresponding to amino acid residues of human NOTCH 2 located near the N-terminal sequence of the cleaved intracellular domain (NICD).
<b>Anti-NOTCH 4</b>	N44-363R	Rabbit, IgG	ELISA, IHC	Rabbit polyclonal NOTCH4 antibody was raised against a peptide from the C terminal residues of human NOTCH4.

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-p16INK4A</b>	C30-363R	Rabbit, IgG	WB, ELISA, IF, IHC	p16INK4A antibody was raised against a 18 amino acid synthetic peptide from near the amino terminus of human p16INK4A
<b>Anti-p16INK4A (CDKN2A)</b>	C30-362R	Rabbit, Antiserum	ELISA, IP, WB	The antibody was produced against human p16 protein (missing first 7 amino acids)
<b>Anti-p21CIP1</b>	C27-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 143-147 of human p21CIP1
<b>Anti-p21CIP1 (p21/WAF1)</b>	C27-61M	Mouse, IgG	WB, ELISA, IP	Produced by repeated immunizations with recombinant human p21 fusion protein.
<b>Anti-p27KIP1 (CDKN1B)</b>	C28-362R	Rabbit, Antiserum	ELISA, IP, WB	The antibody was produced against full length recombinant human p27 protein
<b>Anti-p28ING5</b>	I32-363EG	Goat, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to amino acids 127-140 of human p28 ING5 protein (inhibitor of growth family, member 5)
<b>Anti-p29ING4</b>	I32-363DG	Goat, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to amino acids 164-175 of human p29 ING4 protein (inhibitor of growth family, member 4)
<b>Anti-p33ING1</b>	I32-363G	Goat, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to amino acids 285-296 of Human p33 ING1 protein (inhibitor of growth family, member 1)
<b>Anti-p38 alpha</b>	M39-60M	Mouse, Monoclonal, 9F12	WB	Recombinant full length active p38α protein
<b>Anti-p47ING3</b>	I32-363CG	Goat, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to amino acids 294-304 of Human ING3 protein (Inhibitor of growth family, member 3).
<b>Anti-p53</b>	P05-63R	Rabbit	WB	Human p53 peptide (15-40aa)-KLH conjugates
<b>Anti-p53</b>	P05-60M	Mouse, IgG2a	ELISA, ChIP, IHC, IP, WB	The antibody was produced against recombinant human p53 peptide
<b>Anti-p57KIP2</b>	C29-363R	Rabbit, IgG	WB, ELISA, IF	p57KIP2 antibody was raised against a 17 amino acid synthetic peptide near the amino terminus of human p57KIP2
<b>Anti-p70S6K</b>	R21-63R	Rabbit	WB	C-ISKRPEHLRMNL peptide KLH conjugates
<b>Anti-p70S6K</b>	R21-63BR	Rabbit	WB	Synthetic peptide corresponding to amino acids 409-413 of human p70S6K (protein accession number P23443)
<b>Anti-PACRG</b>	P216-363R	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to amino acids 204-215 of human PACRG protein
<b>Anti-PAK2</b>	P03-63BR	Rabbit, IgG	WB, ELISA, IF, IHC	PAK2 antibody was raised against a 14 amino acid synthetic peptide from near the amino terminus of human PAK2
<b>Anti-PAK2</b>	P03-63R	Rabbit, IgG	WB, ELISA, IF, IHC	PAK2 antibody was raised against a 14 amino acid synthetic peptide from near the carboxy terminus of human PAK2
<b>Anti-PAK4</b>	P05-163R	Rabbit, IgG	WB, ELISA, ICC, IF	PAK4 antibody was raised against a 13 amino acid synthetic peptide from near the center of human PAK4

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-PAK6</b>	P06-63R	Rabbit, IgG	WB, ELISA, IF, IHC	PAK6 antibody was raised against a 13 amino acid synthetic peptide from near the center of human PAK6
<b>Anti-PAK7</b>	P07-63R	Rabbit, IgG	WB, ELISA, ICC, IF	PAK7 antibody was raised against a 13 amino acid synthetic peptide from near the center of human PAK7
<b>Anti-PCNA</b>	P11-363R	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to an internal region of human PCNA
<b>Anti-PDGFB</b>	P842-63R	Rabbit, IgG	WB, ELISA	Rabbit polyclonal PDGFB antibody was raised against a 19 amino acid peptide near the carboxy terminus of human PDGFB.
<b>Anti-PDHK1</b>	P57-63R	Rabbit, IgG	ELISA, IHC, WB	The antibody was produced against synthesized peptide corresponding to a region near the C-Terminal of human PDK-1 protein
<b>Anti-PGCC1</b>	P845-363R	Rabbit, IgG	WB, ELISA	PGCC1 antibody was raised against an 18 amino acid synthetic peptide from near the carboxy terminus of human PGCC1.
<b>Anti-PHLPP1</b>	P71-263R	Rabbit, IgG	WB, ELISA	PHLPP1 antibody was raised against an 18 amino acid synthetic peptide near the amino terminus of human PHLPP1
<b>Anti-PHLPP2</b>	P72-263R	Rabbit, IgG	WB, ELISA	PHLPP2 antibody was raised against a 20 amino acid synthetic peptide near the carboxy terminus of human PHLPP2
<b>Anti-PIAS1</b>	P280-63R	Rabbit, IgG	WB, ELISA, IF, IHC	PIAS1 antibody was raised against a 15 amino acid synthetic peptide near the carboxy terminus of human PIAS1
<b>Anti-PIAS2</b>	P281-63R	Rabbit, IgG	WB, ELISA, IF, IHC	PIAS1 antibody was raised against a 15 amino acid synthetic peptide near the carboxy terminus of human PIAS1
<b>Anti-PIAS3</b>	P282-63R	Rabbit, IgG	WB, ELISA, IF, IHC	PIAS3 antibody was raised against a 12 amino acid synthetic peptide near the carboxy terminus of human PIAS3
<b>Anti-PIAS4</b>	P283-63R	Rabbit, IgG	WB, ELISA, IF, IHC	PIAS4 antibody was raised against a 19 amino acid synthetic peptide near the amino terminus of human PIAS4
<b>Anti-PKAc</b>	P51-63R	Rabbit	WB	Protein Code: P17612
<b>Anti-PKC delta/theta</b>	P64-63DR	Rabbit	WB	Synthesized peptide derived from C-terminus of human PKCθ protein
<b>Anti-PKC zeta</b>	P75-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 408-412 (TSTFC) of human PKCζ
<b>Anti-PLAC1</b>	P38-363R	Rabbit, IgG	WB, ELISA, ICC, IF	PLAC1 antibody was raised against a 15 amino acid synthetic peptide near the carboxy terminus of human PLAC1
<b>Anti-PLAC2</b>	P38-363BR	Rabbit, IgG	WB, ELISA, ICC, IF	PLAC2 antibody was raised against a 17 amino acid synthetic peptide near the center of human PLAC2
<b>Anti-PLAC3</b>	P38-363CR	Rabbit, IgG	WB, ELISA, ICC, IF	PLAC3 antibody was raised against a 15 amino acid synthetic peptide near the amino terminus of human PLAC3
<b>Anti-PLAC4</b>	P38-363DR	Rabbit, IgG	WB, ELISA, ICC, IF	PLAC4 antibody was raised against a 17 amino acid synthetic peptide near the center of human PLAC4

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-PLC gamma 1</b>	P40-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 769-773 of human PLCγ1
<b>Anti-PTEN</b>	P23-263BR	Rabbit, IgG	WB, ELISA, IF	PTEN antibody was raised against a 16 amino acid synthetic peptide from near the carboxy terminus of human PTEN
<b>Anti-PTEN</b>	P23-263R	Rabbit, IgG	WB, ELISA	PTEN antibody was raised against a 16 amino acid synthetic peptide from near the carboxy terminus of human PTEN
<b>Anti-PTK7</b>	P95-63R	Rabbit, IgG	WB, ELISA, IF, IHC	PTK7 antibody was raised against a 18 amino acid synthetic peptide from near the carboxy terminus of human PTK7
<b>Anti-PTK7</b>	P95-63BR	Rabbit, IgG	WB, ELISA, IHC	PTK7 antibody was raised against a 15 amino acid synthetic peptide from near the amino terminus of human PTK7
<b>Anti-PTPN11 (SHP2)</b>	P38-63CR	Rabbit, IgG	WB, ELISA, IF	PTPN11 antibody was raised against a 14 amino acid synthetic peptide from near the carboxy terminus of human PTPN11.
<b>Anti-PTPN11 (SHP2)</b>	P38-63BR	Rabbit, IgG	WB, ELISA	PTPN11 antibody was raised against a 14 amino acid synthetic peptide from near the carboxy terminus of human PTPN11.
<b>Anti-PYK2</b>	P92-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 420-424 of human PYK2
<b>Anti-QIK</b>	S15-63R	Rabbit, IgG	WB, ELISA	QIK antibody was raised against a 16 amino acid synthetic peptide near the center of human QIK
<b>Anti-RAF1</b>	R01-63R	Rabbit, IgG	WB, ELISA	RAF1 antibody was raised against a 19 amino acid synthetic peptide near the amino terminus of human RAF1
<b>Anti-Raptor</b>	R47-363R	Rabbit, IgG	WB, ELISA, ICC	Raptor (CT) polyclonal antibody was raised against a 16 amino acid synthetic peptide from near the carboxy terminus of mouse Raptor
<b>Anti-Raptor</b>	R47-363BR	Rabbit, IgG	WB, ELISA, ICC, IF	Raptor polyclonal antibody was raised against a 13 amino acid synthetic peptide from near the amino-terminus of human Raptor
<b>Anti-RB2 (p130)</b>	R50-363R	Rabbit, Antiserum	ELISA, IHC, IP, WB	The antibody was produced against synthesized peptide corresponding to a region near the C-Term of the human protein conjugated to KLH
<b>Anti-Rel A</b>	R29-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 274-278 of human Rel A
<b>Anti-Rel A</b>	R29-63BR	Rabbit	WB	Synthetic peptide corresponding to amino acids 527-531 of human REL A
<b>Anti-RHEB</b>	R42-363R	Rabbit, IgG	WB, ELISA	Rabbit RHEB polyclonal antibody was raised against a 14 amino acid synthetic peptide from the amino terminus of human RHEB
<b>Anti-RHEB</b>	R42-363BR	Rabbit, IgG	WB, ELISA, IF, IHC	RHEB antibody was raised against a 15 amino acid synthetic peptide from the middle region of human RHEB
<b>Anti-RIPK1</b>	R07-63R	Rabbit, IgG	WB, ELISA, IF, IHC	RIPK1 antibody was raised against a 15 amino acid synthetic peptide from near the amino terminus of human RIPK1

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Anti-RIPK2	R08-63R	Rabbit, IgG	WB, ELISA, ICC	RIPK2 antibody was raised against a peptide corresponding to amino acids near the amino terminus of human RIPK2
Anti-RIPK2	R08-63BR	Rabbit, IgG	WB, ELISA, ICC, IF	RIPK2 antibody was raised against a peptide corresponding to amino acids near the amino terminus of human RIPK2.
Anti-RIPK3	R09-63R	Rabbit, IgG	WB, ELISA, IF, IHC	RIPK3 antibody was raised against a peptide corresponding to amino acids 473 to 486 of murine RIPK3
Anti-ROCK1	R10-63R	Rabbit, IgG	WB, ELISA, ICC, IF	ROCK1 antibody was raised against a 15 amino acid synthetic peptide near the amino terminus of human ROCK1
Anti-ROCK2	R11-63R	Rabbit, IgG	WB, ELISA, IF, IHC	ROCK2 antibody was raised against a 15 amino acid synthetic peptide near the center of human ROCK2
Anti-RSK (pan)	R15-63BR	Rabbit	WB	Synthetic peptide corresponding to amino acids 346-350 of mouse p90 RSK (protein accession number P18653)
Anti-RSK1	R15-63R	Rabbit	WB	C-LAQRVRKLPSTTL peptide KLH conjugates
Anti-RSK1	R15-63CR	Rabbit, IgG	WB, ELISA, ICC, IF	RSK1 antibody was raised against a 16 amino acid synthetic peptide from near the carboxy terminus of human RSK1
Anti-RSK3	R16-63R	Rabbit, IgG	WB	Synthetic peptide corresponding to amino acid residues from the C-terminal region conjugated to KLH
Anti-RSL1D1	R76-363R	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to an internal sequence of human PBK1.
Anti-SGK1	S06-63R	Rabbit, IgG	WB, ELISA, IF, IHC	SGK1 antibody was raised against an 18 amino acid synthetic peptide near the carboxy terminus of human SGK1
Anti-SH2D1A (SAP)	S03-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 113-128 of human SH2D1A (SAP)
Anti-SHP2	P38-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 578-582 of human SHP2
Anti-SIK1	S14-63R	Rabbit, IgG	WB, ELISA, IHC	SIK1 antibody was raised against an 18 amino acid synthetic peptide near the center of human SIK1
Anti-SIK3	S12-63R	Rabbit, IgG	WB, ELISA, IF	SIK3 antibody was raised against an 18 amino acid synthetic peptide near the carboxy terminus of human SIK3
Anti-SIRT1	S35-463R	Rabbit, IgG	WB, ELISA, IF	SIRT1 antibody was raised against a 19 amino acid synthetic peptide near the amino terminus of human SIRT1
Anti-SIRT2	S36-463BR	Rabbit, IgG	WB, ELISA, IF, IHC	SIRT2 antibody was raised against a 17 amino acid synthetic peptide near the amino terminus of the human SIRT2
Anti-SIRT2	S36-463R	Rabbit, IgG	WB, ELISA, IF, IHC	SIRT2 antibody was raised against a 19 amino acid synthetic peptide near the carboxy terminus of the human SIRT2
Anti-SIRT3	S37-463R	Rabbit, IgG	WB, ELISA	SIRT3 antibody was raised against a 17 amino acid synthetic peptide near the center of human SIRT3

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-SIRT4</b>	S38-463R	Rabbit, IgG	WB, ELISA	SIRT4 antibody was raised against a 17 amino acid synthetic peptide near the amino terminus of human SIRT4
<b>Anti-SIRT5</b>	S39-463R	Rabbit, IgG	WB, ELISA	SIRT5 antibody was raised against a 17 amino acid synthetic peptide near the carboxy terminus of human SIRT5
<b>Anti-SIRT6</b>	S40-463R	Rabbit, IgG	WB, ELISA	SIRT6 antibody was raised against a 17 amino acid synthetic peptide near the carboxy terminus of human SIRT6
<b>Anti-SIRT7</b>	S41-463R	Rabbit, IgG	WB, ELISA	SIRT7 antibody was raised against a 19 amino acid synthetic peptide near the amino terminus of human SIRT7
<b>Anti-SMAD1</b>	S10-363R	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to the region of amino acids containing Ser206 of human SAMD1 protein
<b>Anti-SMAD2</b>	S11-363R	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to an internal region of human SMAD2 protein.
<b>Anti-SMC1</b>	S19-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 955-959 of human SMC1 (Q14683)
<b>Anti-SNCA</b>	S22-360M	Mouse, IgG1, 3H9	WB, IF	Full length human alpha synuclein expressed in E. Coli.
<b>Anti-SOX2</b>	S832-361M	Mouse, IgG	WB, ELISA, IHC	Recombinant human SOX2 protein is used to produce this monoclonal antibody
<b>Anti-Src</b>	S19-63BR	Rabbit, IgG	WB, IHC	The antiserum was produced against synthesized non-phosphopeptide derived from human Src around the phosphorylation site of tyrosine 529 (P-Q-YP-Q-P).
<b>Anti-STAT1 alpha</b>	S52-63BR	Rabbit, IgG	WB, ELISA, ICC, IP	STAT1 antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human STAT1 alpha. The sequences differ from the murine corresponding sequences by four amino acids.
<b>Anti-STAT3</b>	S54-63CR	Rabbit, IgG	WB, ELISA	Rabbit polyclonal STAT3 antibody was raised against a 19 amino acid peptide near the amino terminus of human STAT3
<b>Anti-Stathmin</b>	S71-363R	Rabbit, IgG	WB, ELISA, IF, IHC	Stathmin antibody was raised against a 20 amino acid synthetic peptide from near the carboxy terminus of human Stathmin
<b>Anti-Stathmin</b>	S71-363BR	Rabbit, IgG	WB, ELISA, IF, IHC	Stathmin antibody was raised against a 20 amino acid synthetic peptide from near the carboxy terminus of human Stathmin
<b>Anti-STK39 (STLK3)</b>	S51-63R	Rabbit, IgG	WB, ELISA, IF, IHC	STK39 antibody was raised against a 19 amino acid synthetic peptide from near the carboxy terminus of human STK39
<b>Anti-STK39 (STLK3)</b>	S51-63BR	Rabbit, IgG	WB, ELISA, IHC	STK39 antibody was raised against a 18 amino acid synthetic peptide from near the center of human STK39
<b>Anti-SUMO1</b>	S293-63R	Rabbit, IgG	WB, ELISA, IF, IHC	SUMO1 antibody was raised against a 14 amino acid synthetic peptide from near the amino terminus of human SUMO1
<b>Anti-SUMO2/3</b>	S294-63BR	Rabbit, IgG	WB, ELISA, IF, IHC	SUMO2/3 antibody was raised against a 15 amino acid synthetic peptide near the carboxy terminus of human SUMO2.

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Anti-SUMO3	S295-63R	Rabbit, IgG	WB, ELISA, IF, IHC	SUMO3 antibody was raised against a 13 amino acid synthetic peptide near the carboxy terminus of human SUMO3.
Anti-TAB1	T01-363R	Rabbit, IgG	WB, ELISA, ICC, IF	TAB1 antibody was raised against a synthetic peptide corresponding to 13 amino acids in the center of human TAB1
Anti-TAK1	M15-63R	Rabbit, IgG	WB, ELISA, IF, IHC	TAK1 antibody was raised against a synthetic peptide corresponding to 14 amino acids near the amino-terminus of human TAK1 isoform a
Anti-TBK1	T02-63R	Rabbit, IgG	WB, ELISA, ICC	TBK1 antibody was raised against a synthetic peptide corresponding to 17 amino acids form near the carboxy terminus of human TBK1.
Anti-TET3	T638-63R	Rabbit, IgG	WB, ELISA, IF	TET3 antibody was raised against a 17 amino acid synthetic peptide near the carboxy terminus of human TET3
Anti-TRAF2	T57-363BR	Rabbit, IgG	WB, ELISA, IF, IHC	TRAF2 antibody was raised against a 17 amino acid synthetic peptide from near the carboxy terminus of human TRAF2
Anti-TRAF2	T57-363R	Rabbit, IgG	WB, ELISA, IHC	TRAF2 antibody was raised against a 17 amino acid synthetic peptide from near the carboxy terminus of human TRAF2
Anti-TRAF3	T58-363R	Rabbit, IgG	WB, ELISA, IF	TRAF3 antibody was raised against a 16 amino acid synthetic peptide from near the carboxy terminus of human TRAF3
Anti-TRAF3	T58-363BR	Rabbit, IgG	WB, ELISA, ICC	TRAF3 antibody was raised against a 15 amino acid synthetic peptide from near the amino terminus of human TRAF3
Anti-Trail	T880-63R	Rabbit, IgG	WB, ELISA, IHC	TRAIL antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human TRAIL.
Anti-TRKA	N16-63R	Rabbit, IgG	IF	The antibody was produced against synthesized non-phosphopeptide derived from human TRKA around the phosphorylation site of tyrosine791 (P-V-YP-L-D).
Anti-TSC1	T71-363R	Rabbit, IgG	WB, ELISA	Rabbit TSC1 (CT) polyclonal antibody was raised against a 15 amino acid synthetic peptide from the carboxy terminus of human TSC1
Anti-TSC1	T71-363BR	Rabbit, IgG	WB, ELISA, ICC, IF	TSC1 antibody was raised against a 15 amino acid synthetic peptide from the middle region of human TSC1. TSC1 antibody will recognize both isoforms of TSC1
Anti-TSC2	T72-363R	Rabbit, IgG	WB, ELISA	TSC2 antibody was raised against a 15 amino acid synthetic peptide from near the carboxy terminus of human TSC2
Anti-TSC2	T72-363BR	Rabbit, IgG	WB, ELISA	TSC2 (NT) was raised against a 14 amino acid synthetic peptide from near the amino terminus of human TSC2
Anti-TTBK1	T18-63R	Rabbit, IgG	WB, ELISA, IF, IHC	TTBK1 antibody was raised against a 16 amino acid synthetic peptide near the carboxy terminus of human TTBK1
Anti-TTBK2	T18-63BR	Rabbit, IgG	ELISA, IHC	TTBK2 antibody was raised against an 18 amino acid synthetic peptide near the carboxy terminus of human TTBK2

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Anti-Tubulin alpha	T91-63R	Rabbit, IgG	WB, ELISA	Tubulin alpha antibody was raised against a 16 amino acid peptide near the amino terminus of human Tubulin alpha
Anti-Tubulin beta	T92-63R	Rabbit	WB	Synthetic peptide derived from $\beta$ -tubulin of many animal species
Anti-TYK2 (JTK1)	T21-63R	Rabbit, IgG	WB, ELISA	Rabbit polyclonal TYK2 antibody was raised against a 17 amino acid peptide near the amino terminus of human TYK2
Anti-Tyrosine Hydroxylase	H492-363R	Rabbit, IgG	WB, IF, IHC	SDS-denatured rat tyrosine hydroxylase, purified from pheochromocytoma.
Anti-UBA1 (UBE1)	U201-62R	Rabbit, IgG	ELISA, IHC, WB	The antibody was produced against recombinant protein corresponding to full-length human Ubiquitin Activating Enzyme E1
Anti-UBA2 (SAE2)	U202-63R	Rabbit, IgG	WB, ELISA, ICC, IF	UBA2 antibody was raised against an 18 amino acid synthetic peptide near the carboxy terminus of human UBA2
Anti-UBC9	U19-63R	Rabbit	WB	Synthetic peptide of c-terminal amino acid sequence of human UBC9 (P63279)
Anti-UBE2N (UBC13)	U231-63R	Rabbit, IgG	WB, ELISA, IF, IHC	UBE2N (UBC13) antibody was raised against a peptide corresponding to 15 amino acids near the C-terminus of human UBE2N (UBC13).
Anti-UBE2V1 (UEV1A)	U238-63R	Rabbit, IgG	WB, ELISA	UBE2V1 (UEV1A) antibody was raised against a peptide corresponding to 15 amino acids near the C-terminus of human UBE2V1 (UEV1A)
Anti-UBLE1A (SAE1)	U208-63R	Rabbit, IgG	WB, ELISA	UBLE1A antibody was raised against a 20 amino acid synthetic peptide near the carboxy terminus of human UBLE1A
Anti-VASP	V12-63R	Rabbit	WB	Synthetic peptide corresponding to amino acids 155-159 of human VASP
Anti-VAV	V14-63R	Rabbit	WB, IP	Synthetic peptide corresponding to amino acids 172-176 of human VAV
Anti-Vimentin	V49-362C	Chicken, IgY	WB, IF	Recombinant human vimentin purified from E. coli
Anti-VLK	V05-163R	Rabbit, IgG	WB, ELISA, IF, IHC	VLK antibody was raised against a 15 amino acid synthetic peptide from near the center of human VLK
Anti-XIAP	X08-63R	Rabbit, IgG	WB, ELISA, IF, IHC	XIAP antibody was raised against a synthetic peptide corresponding to 13 amino acids at the C-terminus of human XIAP
Anti-ZAP70	Z02-61M	Mouse, IgG	WB	Recombinant Human ZAP70
Anti-ZAP70	Z02-63R	Rabbit, IgG	WB, ELISA	Raised against a recombinant human ZAP70 protein 168-421aa
Anti-ZIP11	Z45-363R	Rabbit, IgG	WB, ELISA, IHC	ZIP11 antibody was raised against a 15 amino acid synthetic peptide near the center of the human ZIP11
Tau Antibody	T08-62BR	Rabbit	WB, IHC, ELISA	Recombinant Human Tau 2NR4 P301S Fibril
Tau Antibody	T08-63R	Rabbit	WB	Recombinant Human Tau AA297-391 (dGAE) Fibril
Tau Antibody	T08-62R	Rabbit	WB, IHC, ELISA	Recombinant Human Tau 2NR4 P301S Fibril

# Primary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Tau Antibody (pSer202/pThr205), AH36</b>	T08-65R	Rabbit / IgG / AH36	WB, ICC/IF, IHC, DB, ELISA	Synthetic peptide of Human Phospho Tau (Ser202/Thr205)
<b>Tau Antibody, 1D5</b>	T08-60M	Mouse / IgG1 / 1D5	WB, DB, ICC/IF, ELISA	Human Recombinant Tau441 (2N4R), P301S mutant Protein Pre-formed fibrils
<b>Tau Antibody, 3D4</b>	T08-60BM	Mouse / IgG1 / 3D4	WB, DB, ICC/IF, ELISA	Human Recombinant Tau441, P301S mutant Protein Pre-formed fibrils

# Secondary Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-GOAT, HRP Conjugated</b>	G34-62DC	Chicken, IgG		The antibody was produced against goat IgG whole molecule
<b>Anti-Mouse HRP</b>	G32-62G	Goat Polyclonal	WB, ELISA	Affinity purified mouse immunoglobulins
<b>Anti-MOUSE, AP Conjugated</b>	M32-62ER	Rabbit, IgG	IHC, ELISA, WB	The antibody was produced against mouse IgG whole molecule
<b>Anti-MOUSE, Biotin Conjugated</b>	M32-62BG	Goat, IgG	WB, ELISA, IHC	The antibody was produced against mouse IgG whole molecule in goat.
<b>Anti-MOUSE, Fluorescein Conjugated</b>	M32-62CG	Goat, IgG	ELISA, IF Microscopy, Flow Cytometry	The antibody was produced against mouse IgG whole molecule.
<b>Anti-MOUSE, Rhodamine Conjugate</b>	M32-62FG	Goat, IgG	ELISA, Flow Cytometry, IF Microscopy	The antibody was produced against mouse IgG whole molecule.
<b>Anti-Rabbit HRP</b>	G33-62G	Goat Polyclonal	WB, ELISA	Affinity purified rabbit immunoglobulins
<b>Anti-RABBIT, AP Conjugated</b>	R33-62EG	Goat, IgG	ELISA, IHC, WB	The antibody was produced against Rabbit IgG whole molecule in goat.
<b>Anti-RABBIT, Biotin Conjugated</b>	R33-62BG	Rabbit, IgG	ELISA, IHC, WB	The antibody was produced against Rabbit IgG whole molecule in goats
<b>Anti-RABBIT, Fluorescein Conjugated</b>	R33-62CG	Goat, IgG	ELISA, Flow Cytometry, IF Microscopy	The antibody was produced against Rabbit IgG whole molecule
<b>Anti-RABBIT, Rhodamine Conjugated</b>	R33-62FG	Goat, IgG	ELISA, FC, IF	The antibody was produced against Rabbit IgG whole molecule.
<b>Anti-RAT, HRP Conjugated</b>	A37-62DG	Goat, IgG	ELISA, IHC, WB	The antibody was produced against rat IgG whole molecule.
<b>Anti-SHEEP, HRP Conjugated</b>	S36-62DD	Donkey, IgG	ELISA, IHC, WB	The antibody was produced against synthesized Sheep IgG whole molecule in donkey.

# Modification-Specific Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-5-Methylated Cytosine (5-methyl cytosine)</b>	C92-67S	Sheep, IgG	WB, IP	5-Methylated Cytosine (5-methyl cytosine)
<b>Anti-Acetylated Lysine</b>	L95-66R	Rabbit	WB, ELISA, PI, IHC, IF	Acetylated lysine-KLH conjugates
<b>Anti-Acetylated Lysine, Biotin Conjugated</b>	L95-66BR	Rabbit	WB, ELISA, IF, IHC, IP	Acetylated lysine-KLH conjugates
<b>Anti-Acetylated Lysine, FITC Conjugated</b>	L95-66CR	Rabbit	For direct immunofluorescence assay	Acetylated lysine-KLH conjugates
<b>Anti-Acetylated Lysine, HRP Conjugated</b>	L95-66DR	Rabbit	WB, ELISA, IF, IHC, IP	Acetylated Lysine-KLH conjugates
<b>Anti-Methylated Arginine (NG, NG-Di-methyl)</b>	A96-67R	Rabbit	WB, ELISA, IF, IHC, IP	NG, NG-di-methylated arginine-KLH conjugates
<b>Anti-Methylated Arginine (NG-Mono-methyl)</b>	A95-67R	Rabbit	WB, ELISA, IF, IHC, IP	NG-mono-methylated arginine-KLH conjugates
<b>Anti-Methylated Lysine (Tri- epsilon-N-methyl)</b>	L95-67R	Rabbit	WB, ELISA, IF, IHC, IP	Methylated lysine-KLH conjugates.
<b>Anti-Methylated Lysine (Tri- epsilon-N-methyl), Biotin Conjugated</b>	L95-67BR	Rabbit	WB, ELISA, IF, IHC	Methylated lysine-KLH conjugates.
<b>Anti-Methylated Lysine (Tri- epsilon-N-methyl), FITC Conjugated</b>	L95-67CR	Rabbit	For direct immunofluorescence assay	Methylated lysine-KLH conjugates.
<b>Anti-Methylated Lysine (Tri- epsilon-N-methyl), HRP Conjugated</b>	L95-67DR	Rabbit	WB, ELISA, IF, IHC, IP	Methylated lysine-KLH conjugates.
<b>Anti-Phosphoserine (Anti-pS)</b>	S95-65R	Rabbit	WB, ELISA, IHC, IP	Phosphoserine peptide-KLH conjugates
<b>Anti-Phosphoserine (Anti-pS), Biotin Conjugated</b>	S95-65BR	Rabbit	WB, ELISA	Phosphoserine peptide-KLH conjugates
<b>Anti-Phosphothreonine (Anti-pT)</b>	T95-65R	Rabbit	WB, ELISA, IHC, IP	Phosphothreonine peptide-KLH conjugates
<b>Anti-Phosphothreonine (Anti-pT), Biotin Conjugated</b>	T95-65BR	Rabbit	WB, ELISA	Phosphothreonine peptide-KLH conjugates
<b>Anti-Phosphotyrosine (Anti-pY)</b>	T96-65R	Rabbit	WB, ELISA, IHC, IP	Phosphotyrosine peptide-KLH conjugates
<b>Anti-Phosphotyrosine (Anti-pY), Biotin Conjugated</b>	T96-65BR	Rabbit	WB, ELISA, IHC, IP	Phosphotyrosine peptide-KLH conjugates

# Phospho-Specific Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Alpha Synuclein Antibody (pSer129), J18</b>	S22-65R	Rabbit / IgG / J18	WB, ELISA, ICC, IHC	Human alpha synuclein AA 124-134: AYEMP-pS-EEGYQ-Cys
<b>Anti-phospho-14-3-3 beta (Ser58)</b>	Y71-365R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser58 conjugated to KLH
<b>Anti-phospho-4EBP1 (Thr45)</b>	E35-365R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr45
<b>Anti-phospho-AKT1/2/3 (Ser473)</b>	A16-65BR	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser473
<b>Anti-phospho-AKT1/2/3 (Ser473)</b>	A16-65BM	Mouse, IgG1	ELISA, IHC, IF, WB	The antibody was produced against synthesized peptide corresponding to residues surrounding Ser 473 of human AKT1 protein.
<b>Anti-phospho-AKT1/2/3 (Thr308)</b>	A16-65R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr308
<b>Anti-phospho-AKT1/2/3 (Thr308)</b>	A16-65M	Mouse, IgG	ELISA, IHC, WB	The antibody was produced against synthesized peptide corresponding to residues surrounding Thr 308 of Human AKT1 protein.
<b>Anti-Phospho-APC1 (Ser377)</b>	A87-365R	Rabbit, IgG	ELISA, IHC, WB	The antibody was produced against synthesized peptide corresponding to amino acids 373-382 of human APC1 protein.
<b>Anti-Phospho-APC3 (Ser427)</b>	A89-365BR	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to amino acids 422-430 of human Cdc27 protein.
<b>Anti-Phospho-APC3 (Thr 244)</b>	A89-365R	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to amino acids 239-240 of Human Cdc27 protein.
<b>Anti-Phospho-ASK1 (Ser83)</b>	M13-65BR	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to amino acids 76-87 on human ASK1 protein.
<b>Anti-phospho-ASK1 (Ser966)</b>	M13-65R	Rabbit, IgG	IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser966
<b>Anti-phospho-ATF2 (Ser112)</b>	A10-565R	Rabbit, IgG	IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser112
<b>Anti-phospho-ATF2 (Ser490 Ser498)</b>	A10-565BR	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser490/498 conjugated to KLH
<b>Anti-Phospho-ATM (Ser1981)</b>	A26-65S	Sheep, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to a region near serine 1981 of human ATM conjugated to KLH.
<b>Anti-Phospho-AURORA B (Thr 232)</b>	A31-65R	Rabbit, IgG	ELISA, IF, WB	The antibody was produced against synthesized peptide corresponding to amino acids 223-234 of Human Aurora Kinase B protein.
<b>Anti-phospho-BRCA1 (Ser1423)</b>	B87-365R	Rabbit, IgG	IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser1423
<b>Anti-phospho-c-KIT (Tyr721)</b>	K06-65R	Rabbit, IgG	WB	The antiserum was produced against synthesized phosphopeptide derived from human c-KIT around the phosphorylation site of tyrosine 721 (N-E-YP-M-D).

# Phospho-Specific Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-phospho-c-KIT (Tyr936)</b>	K06-65BR	Rabbit, IgG	IHC	The antibody was produced against synthesized phosphopeptide derived from human c-KIT around the phosphorylation site of tyrosine 936 (H-I-Y-P-S-N).
<b>Anti-phospho-CAMK2A/2B (Thr286)</b>	C11-65R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr286 conjugated to KLH
<b>Anti-phospho-Catenin beta (Ser33 Ser37)</b>	C06-365R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser33/37 conjugated to KLH
<b>Anti-phospho-CDK1 (Tyr15)</b>	C22-65R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr15 conjugated to KLH
<b>Anti-Phospho-CDK9 (Thr29)</b>	C40-65R	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to residues surrounding Thr29 in the human CDK9 protein.
<b>Anti-phospho-CHK2 (Thr68)</b>	C48-65R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr68
<b>Anti-phospho-CREB1 (Ser133)</b>	C49-565R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser383 conjugated to KLH
<b>Anti-Phospho-Cyclin B1 (Ser126)</b>	C79-365R	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to amino acids 12-131 of Human Cyclin B1 protein
<b>Anti-phospho-EGFR (Thr678)</b>	E10-65BR	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr678
<b>Anti-phospho-EGFR (Thr693)</b>	E10-65CR	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr693
<b>Anti-phospho-EGFR (Tyr1092)</b>	E10-65DR	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr1092
<b>Anti-phospho-EGFR (Tyr1110)</b>	E10-65ER	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr1110
<b>Anti-phospho-EGFR (Tyr1197)</b>	E10-65FR	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr1197
<b>Anti-phospho-EIF2AK2 (Thr446)</b>	P80-65R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr446
<b>Anti-phospho-ELK1 (Ser383)</b>	E64-65R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser383 conjugated to KLH
<b>Anti-phospho-EPHB2 (Tyr469)</b>	E22-65R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr469 conjugated to KLH
<b>Anti-phospho-EPHB2 (Tyr473)</b>	E22-65BR	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr473 conjugated to KLH
<b>Anti-phospho-ERK1/2 (Thr202)</b>	M29-65R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr202
<b>Anti-phospho-ERK1/2 (Thr202/ Tyr204)</b>	M29-652R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr202/Tyr204 conjugated to KLH

# Phospho-Specific Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-phospho-ERK1/2 (Tyr204)</b>	M29-65BR	Rabbit, IgG	WB, IF, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr204
<b>Anti-phospho-GABABR2 (Ser783)</b>	G803-465BR	Rabbit, IgG	WB, IF	Phosphopeptide corresponding to amino acid residues surrounding the phospho-Ser 783 of GABABR2
<b>Anti-phospho-GluR1 (Ser845)</b>	G830-365R	Rabbit, IgG	WB, IHC	Phosphopeptide corresponding to the amino acid residues surrounding the phospho-Ser845 of GluR1.
<b>Anti-Phospho-GSKa (Ser21)</b>	G08-65R	Rabbit, IgG	ELISA, IF, WB	The antibody was produced against synthesized peptide corresponding to amino acids 16-25 of human GSK3 alpha
<b>Anti-phospho-GSK3 beta (Ser9)</b>	G09-652R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser9 conjugated to KLH
<b>Anti-phospho-GSK3 beta (Ser9)</b>	G09-65R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser9
<b>Anti-phospho-HDAC2 (Ser394)</b>	H84-65R	Rabbit, IgG	WB, IF, IHC	Phosphopeptide derived from human HDAC2 around the phosphorylation site of serine 394 (E-D-SP-G-D).
<b>Anti-phospho-HER2 (Tyr877)</b>	E27-65R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr877
<b>Anti-phospho-Histone H2A.X (Ser139)</b>	H11-565R	Rabbit, IgG	WB	The antiserum was produced against synthesized phosphopeptide derived from human Histone H2A.X around the phosphorylation site of serine 139 (Q-A-SP-Q-E).
<b>Anti-phospho-Histone H3.1 (Ser10)</b>	H12-565R	Rabbit, IgG	WB, IHC	The antiserum was produced against synthesized phosphopeptide derived from human Histone H3.1 around the phosphorylation site of serine 10 (R-K-SP-T-G).
<b>Anti-phospho-HTT (Ser421)</b>	H390-65R	Rabbit, IgG	ELISA, IHC, WB	The antibody was produced against synthesized peptide corresponding to amino acids 416-424 of human Huntington protein.
<b>Anti-Phospho-IKKe</b>	I04-65R	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to a region of the protein surrounding phosphorylated Thr501 conjugated to KLH
<b>Anti-Phospho-IRS1 (Ser307)</b>	I40-365CR	Rabbit, IgG	ELISA, IHC, WB	The antibody was produced against synthesized peptide corresponding to amino acids 298-316 of human IRS1 protein.
<b>Anti-phospho-IRS1 (Ser312)</b>	I40-365R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser312
<b>Anti-phospho-IRS1 (Ser639)</b>	I40-365BR	Rabbit, IgG	IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser639
<b>Anti-phospho-JAK1 (Tyr1022)</b>	J01-65R	Rabbit, IgG	IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr1022
<b>Anti-phospho-JAK2 (Tyr1007)</b>	J02-65BR	Rabbit, IgG	IHC	The antiserum was produced against synthesized phosphopeptide derived from human JAK2 around the phosphorylation site of tyrosine 1007 (K-E-YP-Y-K).
<b>Anti-phospho-JAK2 (Tyr221)</b>	J02-65R	Rabbit, IgG	IHC	The antiserum was produced against synthesized phosphopeptide derived from human JAK2 around the phosphorylation site of tyrosine 221 (Q-D-YP-H-I).

# Phospho-Specific Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-phospho-KDR (Tyr1214)</b>	K01-65BR	Rabbit, IgG	IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr1214
<b>Anti-phospho-KDR (Tyr951)</b>	K01-65R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr951
<b>Anti-phospho-MARCKS (Ser152 Ser156)</b>	M02-565R	Rabbit, IgG	WB	corresponding to amino acid residues surrounding Ser152/156 conjugated to KLH
<b>Anti-Phospho-MARK2 (Thr595)</b>	M44-65R	Rabbit, IgG	ELISA, IF, WB	The antibody was produced against synthesized peptide corresponding to amino acids surrounding phosphorylated Thr595 of human MARK2 isoform a.
<b>Anti-Phospho-MDM2 (Ser185)</b>	M45-365R	Rabbit, IgG	ELISA, IP, WB	The antibody was produced against synthesized peptide corresponding to amino acids 177-198 of mouse MDM2
<b>Anti-phospho-MEK1 (Thr292)</b>	M02-65BR	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr292 conjugated to KLH
<b>Anti-phospho-MEK1 (Thr386)</b>	M02-65CR	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr386 conjugated to KLH
<b>Anti-phospho-MEK1/2 (Ser218 Ser222)</b>	M02-652R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser218/222 conjugated to KLH
<b>Anti-phospho-MEK1/2 (Ser222)</b>	M02-65R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser222
<b>Anti-phospho-MEK5 (Ser311 Ser315)</b>	M06-65R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser311/Thr315 conjugated to KLH
<b>Anti-phospho-MET (Tyr1349 Tyr1356)</b>	M52-65R	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to residues surrounding Tyr1349 and Tyr1356 of human c-MET protein.
<b>Anti-phospho-mTOR (Ser2448)</b>	F17-65R	Rabbit, IgG	IHC	The antiserum was produced against synthesized phosphopeptide derived from human mTOR around the phosphorylation site of serine 2448 (T-D-SP-Y-S).
<b>Anti-phospho-p38 alpha (Thr180 Tyr182)</b>	M39-652R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr180/Tyr182 conjugated to KLH
<b>Anti-phospho-p38 alpha (Tyr182)</b>	M39-65R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr182
<b>Anti-phospho-p53 (Ser392)</b>	P05-363R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser392 conjugated to KLH
<b>Anti-phospho-PAK1 (Thr212)</b>	P02-65BR	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr212
<b>Anti-phospho-PAK1/2/3 (Thr402)</b>	P02-652R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr402 conjugated to KLH
<b>Anti-phospho-PAK1/2/3 (Thr423)</b>	P02-653R	Rabbit, IgG	ELISA, WB	The antibody was produced against human PAK1/2/3 phospho peptide corresponding to a region in the protein conjugated to KLH.
<b>Anti-phospho-Parkin (Ser101)</b>	P215-65R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser101 conjugated to KLH

# Phospho-Specific Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-phospho-Parkin (Ser378)</b>	P215-65BR	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser378 conjugated to KLH
<b>Anti-phospho-PDK1 (Ser241)</b>	P14-65R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser241
<b>Anti-phospho-PLCG1 (Tyr783)</b>	P40-465R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr783
<b>Anti-phospho-PLCG2 (Tyr753)</b>	P41-465R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr753
<b>Anti-phospho-PLK1 (Thr210)</b>	P41-65R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr210 conjugated to KLH
<b>Anti-phospho-RAF1 (Ser301)</b>	R01-65R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser301 conjugated to KLH
<b>Anti-phospho-RAF1 (Ser642)</b>	R01-65BR	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser642 conjugated to KLH
<b>Anti-Phospho-Rel A (ser276)</b>	R29-365R	Rabbit, Antiserum	ELISA, IHC, WB	The antibody was produced against synthesized peptide corresponding to a region near phosphorylated Ser276 of the protein conjugated to KLH.
<b>Anti-Phospho-ROCK2 (Tyr256)</b>	R11-65R	Rabbit, IgG	ELISA, IHC, WB	The antibody was produced against synthesized peptide corresponding to residues surrounding Tyrosine 256 of human ROCK-2
<b>Anti-Phospho-RSK1 (Ser732)</b>	R15-65R	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to a region near the C-Terminal of the human RSK1 protein
<b>Anti-Phospho-SMAD1 (Ser206)</b>	S10-365R	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to the region of amino acids containing Ser206 of the SMAD1 protein
<b>Anti-Phospho-SMAD3 (Ser423)</b>	S12-365R	Rabbit, IgG	ELISA, IHC, WB	The antibody was produced against synthesized synthesized peptide corresponding to amino acids 417-425 of human SMAD3 protein.
<b>Anti-Phospho-SMAD3 (Thr179)</b>	S12-365BR	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to an internal region of human SMAD3 protein surrounding Thr179.
<b>Anti-phospho-SNCA (Ser129)</b>	S22-365R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser129 of Rat SNCA conjugated to KLH
<b>Anti-phospho-Src (Tyr529)</b>	S19-65BR	Rabbit, IgG	WB, IHC	The antiserum was produced against synthesized non-phosphopeptide derived from human Src around the phosphorylation site of tyrosine 529 (P-Q-YP-Q-P).
<b>Anti-Phospho-STAT2 (Tyr690)</b>	S53-365R	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to a region near the N-Terminus of human STAT2 protein.
<b>Anti-phospho-STAT3 (Ser727)</b>	S54-365BR	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser727

# Phospho-Specific Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
Anti-phospho-STAT3 (Tyr705)	S54-365R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr705
Anti-Phospho-STAT5 (Tyr694)	S56-365BM	Mouse, IgG1	ELISA, IHC, IF, WB	The antibody was produced against synthesized peptide corresponding to residues surrounding Tyr694 of mouse STAT5a protein.
Anti-phospho-STAT5A (Ser780)	S56-365R	Rabbit, IgG	WB, IHC	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser780
Anti-Phospho-STAT6 (Tyr641)	S57-365R	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to residues surrounding Tyr641 of human STAT6 protein
Anti-phospho-STK10-a (Ser482 Ser486 Ser490)	S29-165R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser482/486/490 conjugated to KLH
Anti-phospho-Tau (Ser214)	T08-365GR	Rabbit, IgG	WB	The antiserum was produced against synthesized phosphopeptide derived from human Tau around the phosphorylation site of serine 235 (P-K-SP-P-S)
Anti-phospho-Tau (Ser235)	T08-365FR	Rabbit, IgG	WB	The antiserum was produced against synthesized phosphopeptide derived from human Tau around the phosphorylation site of serine 235 (P-K-SP-P-S)
Anti-phospho-Tau (Ser262)	T08-365IR	Rabbit, IgG	WB	The antiserum was produced against synthesized phosphopeptide derived from human Tau around the phosphorylation site of serine 262 (I-G-SP-T-E)
Anti-phospho-Tau (Ser356)	T08-365MR	Rabbit, IgG	WB, IF	The antibody was produced against synthesized phosphopeptide derived from human Tau around the phosphorylation site of serine 356 (I-G-SP-L-D)
Anti-phospho-Tau (Ser396)	T08-365ER	Rabbit, IgG	WB	The antiserum was produced against synthesized phosphopeptide derived from human Tau around the phosphorylation site of serine 396 (Y-K-SP-P-V)
Anti-phospho-Tau (Ser404)	T08-365HR	Rabbit, IgG	WB	The antiserum was produced against synthesized phosphopeptide derived from human Tau around the phosphorylation site of serine 404 (D-T-SP-P-R)
Anti-phospho-Tau (Ser416)	T08-365DR	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser416 conjugated to KLH
Anti-phospho-Tau (Ser422)	T08-365KR	Rabbit, IgG	WB	The antibody was produced against synthesized phosphopeptide derived from human Tau around the phosphorylation site of serine 422 (V-D-SP-P-Q)
Anti-phospho-Tau (Thr181)	T08-365R	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr181
Anti-phospho-Tau (Thr205)	T08-365BR	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr205
Anti-phospho-Tau (Thr212)	T08-365JR	Rabbit, IgG	WB	The antibody was produced against synthesized phosphopeptide derived from human Tau around the phosphorylation site of threonine 212 (S-R-TP-P-S)
Anti-phospho-Tau (Thr231)	T08-365CR	Rabbit, IgG	WB	Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr231

# Phospho-Specific Antibodies

Product Name	Cat# No	Host	Cited Application	Immunogen
<b>Anti-phospho-TRKA (Ser791)</b>	N16-65R	Rabbit, IgG	IF	The antibody was produced against synthesized phosphopeptide derived from human TrkA around the phosphorylation site of tyrosine791 (P-V-YP-L-D).
<b>Anti-Phospho-UBLE1A (SAE1) (Ser185)</b>	U208-65R	Rabbit, IgG	ELISA, WB	The antibody was produced against synthesized peptide corresponding to a region surrounding Serine185 of the human SUMO Activating Enzyme E1 protein
<b>Anti-phospho-ZAP70 (Tyr319)</b>	Z02-65R	Rabbit, IgG	WB	The antiserum was produced against synthesized phosphopeptide derived from human ZAP-70 around the phosphorylation site of tyrosine 319 (S-P-YP-S-D).
<b>Anti-Phospho-APC6 (Thr580)</b>	A92-365R	Rabbit, IgG	ELISA, IF, WB	The antibody was produced against synthesized peptide corresponding to amino acids 575-588 of Human APC6 protein.



# Proteins

# Proteins

## Adaptor Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
FADD Protein	F02-30H	Human	HIS	E.coli	Full Length
GAB1 Protein	G03-30H	Human	HIS	Sf9 Cells	Full Length
GRB2 Protein	G45-30G	Human	GST	E.coli	Full Length
GRB7 Protein	G45-30GG	Human	GST	E.coli	Full Length
NCK Protein	N06-30H	Human	HIS	Sf9 Cells	Full Length
NRBP1 Protein	N24-34G	Human	GST	Sf9 Cells	Full Length
TRAF1 Protein	T56-30G	Human	GST	Sf9 Cells	Full Length
TRAF2 Protein	T57-30G	Human	GST	Sf9 Cells	Full Length

## Apoptosis

Product Name	Cat# No	Species	Tag	Expression System	Sequence
ATG13 Protein	A113-30G	Human	GST	E.coli	Full Length
ATG13 Protein	A113-34G	Human	GST	Sf9 Cells	Full length
ATG4B Protein	A104-34H	Human	HIS	E.coli	Full length
BAD Protein	B03-30G	Human	GST	Sf9 Cells	Full Length
BAG1 Protein	B05-31G	Human	GST	E.coli	72-end
BAX Protein	B14-30G	Human	GST	E.coli	Full Length
BCL2 Protein	B25-30G	Human	GST	E.coli	Full Length
BID Protein	B59-30G	Human	GST	Sf9 Cells	Full Length
XIAP Protein	X08-30G	Human	GST	Sf9 Cells	Full Length

## Cell Cycle Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
CDC7 Protein	C26-30G	Human	GST	Sf9 Cells	Full Length
CDK2 Protein	C29-30G	Human	GST	Sf9 Cells	Full Length
CDK5P25 Protein	C34-31G	Human	GST	Sf9 Cells	98-end
CDK5P35 Protein	C34-30G	Human	GST	Sf9 Cells	Full Length
CDKN3 Protein	C01-34G	Human	GST	Sf9 Cells	Full Length
Cyclin E2 Protein	C82-30BG	Human	GST	Sf9 Cells	Full Length
Cyclin T2 Protein	C86-30BG	Human	GST	Sf9 Cells	Full-Length
Cyclin Y Protein	C87-30G	Human	GST	Sf9 Cells	2-end
CyclinA2 Protein	C78-30BG	Human	GST	Sf9 Cells	Full Length
CyclinC Protein	C80-30G	Human	GST	Sf9 Cells	Full Length
CyclinD1 Protein	C81-30G	Human	GST	E.coli	Full Length
CyclinD1 Protein	C81-30BG	Human	GST	Sf9 Cells	Full Length
CyclinD2 Protein	C81-30B2G	Human	GST	Sf9 Cells	Full length
CyclinD3 Protein	C81-30B3G	Human	GST	Sf9 Cells	Full length
CyclinE1 Protein	C82-30G	Human	GST	Sf9 Cells	Full Length

# Cell Cycle Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
CyclinH Protein	C83-30H	Human	HIS	Sf9 Cells	Full Length
CyclinK Protein	C84-30G	Human	GST	Sf9 Cells	Full Length
CyclinM Protein	C89-30G	Human	GST	Sf9 Cells	Full Length
CyclinO Protein	C85-30G	Human	GST	Sf9 Cells	Full Length
MDM2 (1-118) Protein	M45-31BH	Human	HIS	E.coli	1-118
MDM2 Protein	M45-31G	Human	GST	E.coli	202-end
MDM4 (1-137) Protein	M47-31BH	Human	HIS	E.coli	1-137
p15INK4B Protein	C31-30G	Human	GST	E.coli	Full Length
p16INK4A Protein	C30-30G	Human	GST	E.coli	Full Length
p18INK4C Protein	C32-30G	Human	GST	E.coli	Full Length
p19INK4D Protein	C33-30G	Human	GST	E.coli	Full Length
p21CIP1 Protein	C27-30G	Human	GST	E.coli	Full Length
p27KIP1 Protein	C28-30G	Human	GST	E.coli	Full Length
p33ING1 Protein	I32-30H	Human	HIS	Sf9 Cells	Full Length
p33ING2 Protein	I32-30BG	Human	GST	Sf9 Cells	Full Length
p53 Protein	P05-30G	Human	GST	E.coli	Full Length
p53 Protein	P05-30BG	Human	GST	Hi5 insect cells using baculovirus	Full Length
p63 Protein	P06-30G	Human	GST	Sf9 Cells	Full Length
p73 alpha Protein	P08-30G	Human	GST	Sf9 Cells	Full Length
p73 beta Protein	P08-30BG	Human	GST	Sf9 Cells	Full Length
p73 gamma Protein	P08-30CG	Human	GST	Sf9 Cells	Full Length
PCNA Protein	P11-30H	Human	HIS	E.coli	Full Length

# Cell Stress and Chaperone Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
14-3-3 alpha/beta Protein	Y71-30G	Human	GST	E.coli	Full Length
14-3-3 alpha/beta Protein	Y71-30N	Human	TAG-Free	E.coli	Full Length
14-3-3 alpha/beta Protein	Y71-30H	Human	HIS	Sf9 Cells	Full Length
14-3-3 alpha/beta Protein	Y71-30BH	Mouse	HIS	Sf9 Cells	Full Length
14-3-3 epsilon Protein	Y75-30G	Human	GST	E.coli	Full Length
14-3-3 epsilon Protein	Y75-30N	Human	TAG-Free	E.coli	Full Length
14-3-3 epsilon Protein	Y75-30H	Human	HIS	Sf9 Cells	YWHAE, MDS, MDCR, KCIP-1, 14-3-3E, FLJ45465
14-3-3 epsilon Protein	Y75-30BH	Mouse	HIS	Sf9 Cells	
14-3-3 sigma Protein	Y86-30G	Human	GST	E.coli	Full Length
14-3-3 sigma Protein	Y86-30N	Human	TAG-Free	E.coli	Full Length
14-3-3 sigma Protein	Y86-30H	Human	HIS	Sf9 Cells	Full Length
14-3-3 theta Protein	Y84-30G	Human	GST	E.coli	Full Length
14-3-3 theta Protein	Y84-30N	Human	TAG-Free	E.coli	Full Length
14-3-3 theta Protein	Y84-30H	Human	HIS	Sf9 Cells	Full Length

# Cell Stress and Chaperone Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
14-3-3 theta Protein	Y84-30BH	Mouse	HIS	Sf9 Cells	Full Length
14-3-3 zeta Protein	Y92-30G	Human	GST	E.coli	Full Length
14-3-3 zeta Protein	Y92-30N	Human	TAG-Free	E.coli	Full Length
14-3-3 zeta Protein	Y92-30H	Human	HIS	Sf9 Cells	Full Length
14-3-3 zeta Protein	Y92-30BH	Mouse	HIS	Sf9 Cells	Full Length
CRYAA Protein	C256-30H	Human	HIS	E.coli	Full Length
GST Protein	G52-30U	Artificial	TAG-Free	E.coli	1-226
HO1 Protein	H20-30G	Human	GST	E.coli	Full Length
HO2 Protein	H21-30G	Human	GST	E.coli	Full Length
HSP10 Protein	H29-54G	Human	GST	E.coli	Full Length
HSP27 Protein	H30-54G	Human	GST	E.coli	Full Length
HSP40 Protein	H32-54G	Human	GST	E.coli	Full Length
HSP60 Protein	H33-54G	Human	GST	E.coli	Full Length
HSP70 Protein	H34-54G	Human	GST	E.coli	Full Length
HSP70 Protein	H34-54H	Human	HIS	Sf9 Cells	Full Length
HSP90 alpha Protein	H36-50H	Human	HIS	Sf9 Cells	Full Length
HSP90 beta Protein	H36-54H	Human	HIS	Sf9 Cells	Full Length
Mortalin Protein	M54-31G	Human	GST	Sf9 Cells	47-end
PARK7 (DJ-1) Protein	P219-31H	Human	HIS	E.coli	19-end

# Fluorescent and Color Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
BFP Protein	B620-30H	Bacteria	HIS	E.coli	228-end
eRFP Protein	R620-30BH		HIS	E.coli	Full Length
GFP Protein	E620-30H		HIS	E.coli	Full Length
Red Protein	R620-30CH		HIS	E.coli	Full Length
RFP Protein	R620-30H		HIS	E.coli	Full Length
Y1BFP Protein	B620-30BH		HIS	E.coli	Full Length

# Microtubule and Actin-Associated Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
Alpha Synuclein (A53T) Pre-Formed Fibrils (Type 1)	S22-361N	Human	TAG-Free	E.coli	Full Length
Alpha Synuclein (A53T) Protein	S22-36N	Human	TAG-Free	E.coli	Full Length
Alpha Synuclein Pre-Formed Fibril (Type 1)	S22-381N	Human	TAG-Free	E.coli	Full Length
Alpha Synuclein Pre-Formed Fibril (Type 1)	S22M-381N	Mouse	TAG-Free	E.coli	Full Length
Alpha synuclein Pre-Formed Fibril (Type 2)	S22-382N	Human	TAG-Free	E.coli	Full Length
Alpha Synuclein Protein Monomer	S22M-37N	Mouse	TAG-Free	E.coli	Full Length
Alpha Synuclein Protein Monomer (Type 1)	S22-371N	Human	TAG-Free	E.coli	Full Length
Alpha-synuclein Protein	S22-30H	Human	HIS	E.coli	Full Length

# Microtubule and Actin-Associated Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
Alpha-synuclein Protein Monomer (Type 2)	S22-372N	Human	TAG-Free	E.coli	Full length
Alpha-synuclein, PLK2-phosphorylated	S22-54H	Human	HIS	E.coli	Full Length
Beta-synuclein Protein	S22-30BH	Human	HIS	E.coli	Full Length
Catenin alpha Protein	C05-30G	Human	GST	Sf9 Cells	Full Length
Cofilin 1 Protein	C37-54H	Human	HIS	E.coli	Full Length
Gamma-synuclein Protein	S22-30CH	Human	HIS	E.coli	Full Length
INCENP Protein	I30-31H	Human	HIS	Sf9 Cells	783-end
Stathmin Protein	S71-30H	Human	HIS	E.coli	Full Length
TPM3 Protein	T36-30H	Human	HIS	E.coli	Full Length
TPX2 Protein	T40-30H	Human	HIS	Sf9 Cells	Full Length
Troponin1 Protein	T68-30H	Human	HIS	E.coli	Full Length
Tubulin Beta (TUBB1) Protein	T92-50G	Human	GST	E.coli	Full Length
Vimentin Protein	V49-30H	Human	HIS	E.coli	Full Length

## Regulatory Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
4EBP1 Protein	E35-54H	Human	HIS	Sf9 Cells	Full Length
4EBP1 Protein	E35-54G	Human	GST	E.coli	Full Length
4EBP1 Protein	E35-54BH	Human	HIS	E.coli	Full Length
AMPKG1 Protein	A60-30G	Human	GST	Sf9 Cells	Full Length
AMPKG3 Protein	A60-30CG	Human	GST/ HIS	Sf9 Cells	Full Length
Annexin 5 Protein	A61-30H	Human	HIS	E.coli	Full Length
ATF2 Protein	A10-55G	Human	GST	E.coli	1-254
AXIN1 Protein	A71-30G	Human	GST	Sf9 Cells	Full Length
AXIN2 Protein	A72-30G	Human	GST	Sf9 Cells	Full Length
BRD3 Protein	B17-30G	Human	GST	Sf9 Cells	Full Length
Calmodulin 1 Protein	C01-30H	Human	HIS	E.coli	Full Length
CD19 Protein	C19-35H	Human	HIS	CHO cells	20-291
CRKL Protein	C53-30H	Human	HIS	Sf9 Cells	Full Length
DPYSL2 Protein	D470-30BG	Human	GST	Sf9 Cells	Full Length
EIF2S1 Protein	E23-54H	Human	HIS	E.coli	Full Length
EIF4E Protein	E34-30G	Human	GST	E.coli	Full Length
ENPP2 Protein	E22-35G	Human	GST	Sf9 Cells	36-end
ERF1 Protein	E380-80G	Human	GST	Sf9 Cells	2-end
FKBP1A Protein	F37-30G	Human	GST	E.coli	Full Length
FRS3 Protein	F80-30G	Human	GST	Sf9 Cells	Full Length
GAP43 Protein	G15-30H	Human	HIS	E.coli	Full Length
HSD17B10 Protein	H23-30G	Human	GST	Sf9 Cells	Full Length
HtrA (S234A) Protein	H531-36CH	Streptococcus	HIS	E.coli	33-393
HtrA Protein	H531-30CH	Streptococcus	HIS	E.coli	33-393
HTRA1 (161-379) Protein	H531-32H	Human	HIS	E.coli	161-379
HTRA1 (161-480) Protein	H531-31H	Human	HIS	E.coli	161-480

# Regulatory Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
HTRA1 (S328A) Protein	H531-32BH	Human	HIS	E.coli	161-379
HTRA1 Protein	H531-30H	Human	HIS	Sf9 Cells	23-480
HTRA2 Protein	H532-31H	Human	HIS	E.coli	133-358
HTRA3 Protein	H533-31H	Human	HIS	E.coli	132-353
HTRA4 Protein	H534-31H	Human	HIS	E.coli	159-376
HTRA4 Protein	H534-30H	Human	HIS	Sf9 Cells	32-end
ILK Protein	I06-34G	Human	GST	Sf9 Cells	Full length
IRS1 (C-Term) Protein	I40-31G	Human	GST	Sf9 Cells	600-1245
IRS1 (N-Term) Protein	I40-35G	Human	GST	E.coli	1-355
MBP Protein	M42-54G	Human	GST	E.coli	Full Length
MO25 alpha Protein	M51-30G	Human	GST	Sf9 Cells	Full Length
MO25α Protein	M51-30H	Human	HIS	Sf9 Cells	Full Length
NPM1 Protein	N50-30H	Human	HIS	E.coli	Full Length
NPR1 Protein	N22-35G	Human	GST	Sf9 Cells	495-end
NPR2 Protein	N23-35G	Human	GST	Sf9 Cells	479-end
PI3K (p55 gamma) Protein	P31-30CH	Human	HIS	Sf9 Cells	Full Length
PI3K (p65 alpha) Protein	P31-31H	Human	HIS	Sf9 Cells	1-571
PI3K (p85 alpha) Protein	P31-30H	Human	HIS	Sf9 Cells	Full Length
PI3K (p85 beta) Protein	P31-30BH	Human	HIS	Sf9 Cells	Full Length
POR Protein	P242-30G	Human	GST	Sf9 Cells	Full Length
PP2B Protein	P18-34G	Human	GST	Sf9 Cells	Full Length
PPP2R1A Protein	P162-30G	Human	GST	Sf9 Cells	Full Length
PPP2R1A Protein	P162-30H	Human	HIS	Sf9 Cells	Full Length
PPP2R1B Protein	P162-30BG	Human	GST	Sf9 Cells	Full Length
PRKAR1A Protein	P57-30H	Human	HIS	Sf9 Cells	Full Length
PRKAR1B Protein	P59-30H	Human	HIS	Sf9 Cells	Full Length
PRPF4B Protein	P87-35G	Human	GST	Sf9 Cells	551-end
RanGAP1 Protein	R298-31H	Human	HIS	E.coli	418-end
Rb (C-Term) Protein	R05-55G	Human	GST	E.coli	773-928
Rb (N-Term) Protein	R05-56G	Human	GST	E.coli	48-378
RBBP4 Protein	R314-30G	Human	GST	Sf9 Cells	Full Length
SCYL2 Protein	S03-35G	Human	GST	Sf9 Cells	1-520
SCYL3 Protein	S09-30G	Human	GST	Sf9 Cells	2-end
VAV1 Protein	V14-31G	Human	GST	Sf9 Cells	73-end
WDR48 Protein	W298-31G	Human	GST	Sf9 Cells	2-580

# Transcription Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
ATF1 Protein	A09-54G	Human	GST	Sf9 Cells	Full Length
c-JUN Protein	J05-55G	Human	GST	E.coli	1-79
Catenin beta Protein	C06-30G	Human	GST	Sf9 Cells	Full Length
Catenin beta Protein	C06-30H	Human	HIS	Sf9 Cells	Full length
CBP Protein	C07-31G	Human	GST	E.coli	1319-1710
CREB1 Protein	C49-55G	Rat	GST	E.coli	1-280
ELK1 Protein	E64-30G	Human	GST	Sf9 Cells	Full Length
FOS Protein	F66-30G	Human	GST	Sf9 Cells	Full Length
FOXL2 Protein	F71-30BG	Human	GST	Sf9 Cells	Full Length
HIF1B Protein	H07-30BG	Human	GST	Sf9 Cells	Full Length
HSF1 Protein	H25-30G	Human	GST	E.coli	Full Length
IkBA Protein	I20-30G	Human	GST	E.coli	Full Length
IkBA Protein	I20-31G	Human	GST	E.coli	1-175
IkBB Protein	I21-30G	Human	GST	E.coli	Full Length
MEF2A Protein	M20-34G	Human	GST	Sf9 Cells	Full-length
MYC Protein	M86-30G	Human	GST	Sf9 Cells	Full-length
MYCN Protein	M86-34G	Human	GST	Sf9 Cells	Full Length
NFATC1 Protein	N12-30G	Human	GST	Sf9 Cells	Full Length
NFKB2 Protein	N13-31G	Human	GST	Sf9 Cells	1-454
P300 Protein	P07-31G	Human	GST	Sf9 Cells	1283-1673
Rel B Protein	R30-30H	Human	HIS	Sf9 Cells	Full Length
SCYL1 Protein	S04-35G	Human	GST	Sf9 Cells	1-556
SMAD1 Protein	S10-30G	Human	GST	E.coli	Full Length
SMAD2 (del SXS) Protein	S11-31G	Human	GST	E.coli	C-terminal SXS deletion
SMAD2 Protein	S11-30G	Human	GST	E.coli	Full Length
SMAD3 (del SXS) Protein	S12-31G	Human	GST	E.coli	C-terminal SXS deletion
SMAD3 Protein	S12-30G	Human	GST	E.coli	Full Length
SMAD4 Protein	S13-30G	Human	GST	E.coli	Full Length
SMAD5 Protein	S14-30G	Human	GST	E.coli	Full Length
SMAD9 Protein	S17-30G	Human	GST	E.coli	Full Length
STAT1 alpha Protein	S52-50G	Human	GST	Sf9 Cells	Full Length
STAT1 beta Protein	S52-54G	Human	GST	Sf9 Cells	Full Length
STAT2 Protein	S53-54G	Human	GST	Sf9 Cells	Full Length
STAT3 Protein	S54-54G	Human	GST	Sf9 Cells	Full Length
STAT3 Protein	S54-54BH	Human	HIS	Sf9 Cells	Full Length
STAT4 Protein	S55-54H	Human	HIS	Sf9 Cells	Full Length
STAT5 Protein	S56-54H	Human	HIS	Sf9 Cells	Full Length
STAT5B Protein	S56-54BG	Human	GST	Sf9 Cells	Full Length
STAT5B Protein	S56-54BH	Human	HIS	Sf9 Cells	Full Length
STAT6 Protein	S57-30G	Human	GST	Sf9 Cells	Full Length
TAF1 (C-term) Protein	T28-35BG	Human	GST	Sf9 Cells	1032-end
TAF1 (N-term) Protein	T28-35G	Human	GST	Sf9 Cells	1-561
TDP43 Protein	T74-34G	Human	GST	Sf9 Cells	Full Length

# Tau Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
<b>Tau K18 (P61L) Pre-Formed Fibril</b>	T18-58GN		TAG-Free	E.coli	
<b>Tau K18 (P61L) Protein Monomer</b>	T18-57GN	Human	TAG-Free	E.coli	
<b>Tau Protein Marker</b>	T08-07N				
<b>Tau-316 Protein</b>	T02-54N	Human	Tag-free	E.coli	Full Length
<b>Tau-352 Protein</b>	T03-54N	Human	TAG-Free	E.coli	Full Length
<b>Tau-361 Protein</b>	T09B-54N	Drosophila melano-gaster	Tag-free	E.coli	Full Length
<b>Tau-381 Protein</b>	T04-54N	Human	TAG-Free	E.coli	Full Length
<b>Tau-383 (N352H) Protein</b>	T05-56N	Human	TAG-Free	E.coli	Full Length (N352H)
<b>Tau-383 Protein</b>	T05-54N	Human	TAG-Free	E.coli	Full Length
<b>Tau-410 Protein</b>	T06-54N	Human	TAG-Free	E.coli	Full Length
<b>Tau-412 Protein</b>	T07-54N	Human	TAG-Free	E.coli	Full Length
<b>Tau-412, GSK3beta-phosphorylated</b>	T07-50CN	Human		E.coli	Full length
<b>Tau-441 (1-391) Protein</b>	T08-55CN	Human	TAG-Free	E.coli	1-391
<b>Tau-441 (1-421) Protein</b>	T08-55BN	Human	TAG-Free	E.coli	1-421
<b>Tau-441 (127-421) Protein</b>	T08-55KN	Human		E.coli	127-421
<b>Tau-441 (151-391) Protein</b>	T08-55FN	Human	TAG-Free	E.coli	151-391
<b>Tau-441 (151-421) Protein</b>	T08-55EN	Human	TAG-Free	E.coli	151-421
<b>Tau-441 (151-441) Protein</b>	T08-55DN	Human	TAG-Free	E.coli	151-441
<b>Tau-441 (216-391) Protein</b>	T08-55LN	Human	TAG-Free	E.coli	216-391
<b>Tau-441 (216-391), Biotinylated</b>	T08-55LNB	Human	AVI	E.coli	(216-391)
<b>Tau-441 (231-391) Protein</b>	T08-55IN	Human	TAG-Free	E.coli	231-391
<b>Tau-441 (231-421) Protein</b>	T08-55HN	Human	TAG-Free	E.coli	231-421
<b>Tau-441 (231-441) Protein</b>	T08-55GN	Human	TAG-Free	E.coli	231-441
<b>Tau-441 (244-372) Protein</b>	T08-55N	Human	TAG-Free	E.coli	244-372
<b>Tau-441 (244-372) Protein</b>	T08-55H	Human	HIS	E.coli	244-372
<b>Tau-441 (244-372), Biotinylated</b>	T08-55NB	Human	TAG-Free	E.coli	244-372
<b>Tau-441 (50-391) Protein</b>	T08-55ON	Human	TAG-Free	E.coli	50-391
<b>Tau-441 (50-421) Protein</b>	T08-55NN	Human	TAG-Free	E.coli	50-421
<b>Tau-441 (50-441) Protein</b>	T08-55MN	Human	TAG-Free	E.coli	50-441
<b>Tau-441 (99-441) Protein</b>	T08-55JN	Human	TAG-Free	E.coli	99-441
<b>Tau-441 (A152T) Protein</b>	T08-56VN	Human	TAG-Free	E.coli	A152T
<b>Tau-441 (dK280) Protein</b>	T08-52N	Human	TAG-Free	E.coli	deleted K280
<b>Tau-441 (dN296) Protein</b>	T08-52BN	Human	TAG-Free	E.coli	deleted N296
<b>Tau-441 (G272V) Protein</b>	T08-56DN	Human	TAG-Free	E.coli	G272V
<b>Tau-441 (K257T) Protein</b>	T08-56N	Human	TAG-Free	E.coli	K257T
<b>Tau-441 (L266V) Protein</b>	T08-56CN	Human	TAG-Free	E.coli	L266V
<b>Tau-441 (N279K) Protein</b>	T08-56EN	Human	TAG-Free	E.coli	N279K
<b>Tau-441 (P301L) Protein</b>	T08-56FN	Human	TAG-Free	E.coli	P301L
<b>Tau-441 (P301L), Biotinylated</b>	T08-56FBN	Human		E.coli	Full length
<b>Tau-441 (P301S) Pre-Formed Fibril</b>	T08-58GN	Human	TAG-Free	E.coli	Full Length
<b>Tau-441 (P301S) Protein</b>	T08-56GN	Human	TAG-Free	E.coli	P301S
<b>Tau-441 (P301S) Protein Monomer</b>	T08-57GN	Human	TAG-Free	E.coli	Full legnth
<b>Tau-441 (R406W) Protein</b>	T08-56MN	Human	TAG-Free	E.coli	R406W

# Tau Proteins

Product Name	Cat# No	Species	Tag	Expression System	Sequence
Tau-441 (S198A) Protein	T08-53BN	Human	TAG-Free	E.coli	S198A
Tau-441 (S198E) Protein	T08-53N	Human	TAG-Free	E.coli	S198E
Tau-441 (S199E) Protein	T08-53CN	Human	TAG-Free	E.coli	S199E
Tau-441 (S214A) Protein	T08-53DN	Human	TAG-Free	E.coli	S214A
Tau-441 (S305N) Protein	T08-56HN	Human	TAG-Free	E.coli	S305N
Tau-441 (S352L) Protein	T08-56KN	Human	TAG-Free	E.coli	S352L
Tau-441 (S404A) Protein	T08-53HN	Human	TAG-Free	E.coli	S404A
Tau-441 (S404E) Protein	T08-53GN	Human	TAG-Free	E.coli	S404E
Tau-441 (T217A) Protein	T08-56ON	Human	TAG-Free	E.coli	Full length
Tau-441 (V337M) Protein	T08-56JN	Human	TAG-Free	E.coli	V337M
Tau-441 Protein	T08-54N	Human	TAG-Free	E.coli	Full Length
Tau-441 Protein	T08-54H	Human	HIS	E.coli	Full Length
Tau-441, Biotinylated	T08-54BN	Human	TAG-Free	E.coli	Full Length
Tau-441, Biotinylated	T08-54B2N	Human		E.coli	Full length
Tau-441, BRSK1-phosphorylated	T08-50N	Human	TAG-Free	E.coli	1-441
Tau-441, BRSK2-phosphorylated	T08-50BN	Human	TAG-Free	E.coli	1-441
Tau-441, CAMK2-phosphorylated	T08-50CN	Human	TAG-Free	E.coli	1-441
Tau-441, DYRK1A-phosphorylated	T08-50RN	Human	Tag-free	E.coli	Full Length
Tau-441, DYRK1A-phosphorylated & Biotinylated	T08-50RNB	Human	TAG-Free	E.coli	Full Length
Tau-441, GSK3beta-phosphorylated	T08-50FN	Human	TAG-Free	E.coli	1-441
Tau-441, PHKG2-phosphorylated	T08-50KN	Human	TAG-Free	E.coli	1-441
Tau-441, PKA-phosphorylated	T08-50LN	Human	TAG-Free	E.coli	1-441
Tau-441, TTBK1-phosphorylated	T08-50ON	Human	TAG-Free	E.coli	1-441

# Growth Factors

Product Name	Cat# No	Species	Tag	Expression System	Sequence
4-1BB Receptor (TNFRSF9)	F802-40N	Human		E.coli	
Adiponectin (Acrp30)	A801-40N	Human		E.coli	
AITRL (TNFSF18)	A828-40N	Human		E.coli	
AREG	A875-40N	Human		E.coli	
B-cell Maturation Protein (TNFRSF17)	B819-40N	Human		E.coli	
BAFF-R (TNFRSF13C)	B806-40N	Human		E.coli	
Beta-Defensin 1 (DEFB1)	B823-40N	Human		E.coli	
Beta-Defensin 2 (DEFB4A)	D822-40N	Human		E.coli	
Beta-Defensin 3 (DEFB103A)	D823-40N	Human		E.coli	
Cardiotrophin-1 (CT-1)	C801-40N	Human		E.coli	
CCL2	M821-40N	Human		E.coli	
CCL21	E882-40N	Human		E.coli	
CCL22	M832-40N	Human		E.coli	
CCL3	M841-40N	Human		E.coli	
CCL4	M841-40BN	Human		E.coli	
CD40 Ligand (CD40LG)	C810-40N	Human		E.coli	

# Growth Factors

Product Name	Cat# No	Species	Tag	Expression System	Sequence
CNTF	C851-40N	Human		E.coli	
CXCL1	G871-40N	Human		E.coli	
CXCL11	I822-40N	Human		E.coli	
CXCL12 (SDF-1 alpha)	S821-402N	Human		E.coli	
CXCL12 (SDF-1 beta)	S821-40BN	Human		E.coli	
CXCL14 (BRAK)	M842-40N	Human		E.coli	
CXCL16	C890-40N	Human		E.coli	
CXCL2	G872-40N	Human		E.coli	
DHH	D834-40N	Human		E.coli	
EGF	E801-40N	Human		E.coli	
Eotaxin (CCL11)	E852-40N	Human		E.coli	
EPO	E860-40N	Human		CHO cells	
FGF1	F819-40N	Human		E.coli	
FGF10	F830-40N	Human		E.coli	
FGF16	F836-40N	Human		E.coli	
FGF17	F837-40N	Human		E.coli	
FGF18	F838-40N	Human		E.coli	
FGF2	F820-40N	Human		E.coli	
FGF20	F840-40N	Human		E.coli	
FGF21	F841-40N	Human		E.coli	
FGF3	F823-40N	Human		E.coli	
FGF4	F824-40N	Human		E.coli	
FGF8	F828-40N	Human		E.coli	
Fractalkine (CX3CL1)	F870-40N	Human		E.coli	
G-CSF (CSF3)	G805-40N	Human		E.coli	
Galectin-1 (LGALS1)	G802-40N	Human		E.coli	
GNDF	G831-40N	Human		E.coli	
GH1	G807-402N	Human		E.coli	
GM-CSF (CSF2)	G806-40N	Human		E.coli	
GMFB	G842-40N	Human		E.coli	
GRO gamma (CXCL3)	G873-40N	Human		E.coli	
HB-EGF (Heparin Binding EGF)	H805-40N	Human		E.coli	
HCC-1 (CCL14)	H811-40N	Human		E.coli	
IFNA1 (IFN-alpha1)	I834-40N	Human		E.coli	
IFNA4	I836-40N	Human		E.coli	
IGF1	I811-40N	Human		E.coli	
IGF2	I812-40N	Human		E.coli	
IGFBP3	I813-40N	Human		E.coli	
IL11	I871-40N	Human		E.coli	
IL13	I873-40N	Human		E.coli	
IL15	I875-40N	Human		E.coli	
IL16	I876-40N	Human		E.coli	
IL17A	I877-40N	Human		E.coli	
IL17D	I877-40DN	Human		E.coli	
IL17F	I877-40FN	Human		E.coli	

# Growth Factors

Product Name	Cat# No	Species	Tag	Expression System	Sequence
IL19	I879-40N	Human		E.coli	
IL1A	I861-40N	Human		E.coli	
IL1B	I861-40BN	Human		E.coli	
IL1RN	I860-40N	Human		E.coli	
IL2	I862-40BN	Human		E.coli	
IL22	I882-40N	Human		E.coli	
IL28A	I888-40N	Human		E.coli	
IL3	I863-40N	Human		E.coli	
IL33	I893-40N	Human		E.coli	
IL36A	I896-40N	Human		E.coli	
IL36B	I896-40BN	Human		E.coli	
IL36G	I896-40CN	Human		E.coli	
IL37/IL23	I883-40N	Human		E.coli	
IL4	I864-40N	Human		E.coli	
IL5	I865-40N	Human		E.coli	
IL6	I866-40N	Human		E.coli	
IL7	I867-40N	Human		E.coli	
IL8 (72aa)	I868-40BN	Human		E.coli	
IL8 (77aa)	I868-40CN	Human		E.coli	
IL9	I869-40N	Human		E.coli	
IP-10 (CXCL10)	I838-40BN	Human		E.coli	
KGF (FGF7)	K818-40N	Human		E.coli	
Leptin (LEP)	L833-40BN	Human		E.coli	
LIF	L845-40N	Human		E.coli	
LTA	T861-402N	Human		E.coli	
MANF	M805-40N	Human		E.coli	
Maspin (SerpB5)	M806-40N	Human		E.coli	
MCP-3 (CCL7)	M823-40N	Human		E.coli	
MCP-4 (CCL13)	M824-40N	Human		E.coli	
MIA	M837-40N	Human		E.coli	
Midkine (NEGF2)	M838-40N	Human		E.coli	
MIF	M840-40N	Human		E.coli	
MIP-3 (CCL23)	M843-40N	Human		E.coli	
MIP-4 (CCL18)	M844-40N	Human		E.coli	
MMP2	M846-40N	Human		E.coli	
MSTN	G818-40N	Human		E.coli	
Nodal	N822-40N	Human		E.coli	
OSM	O841-40N	Human		E.coli	
PDGFA	P841-40BN	Human		E.coli	
PDGFB	P842-40N	Human		E.coli	
Persephin (PSPN)	P870-40N	Human		E.coli	
PIGF	P829-40N	Human		E.coli	
Prolactin (PRL)	P852-40N	Human		E.coli	
Resistin (RETN)	R821-40N	Human		E.coli	
S100A4	S801-40N	Human		E.coli	

# Growth Factors

Product Name	Cat# No	Species	Tag	Expression System	Sequence
SERPINF1	P843-40N	Human		E.coli	
SHH	S851-40N	Human		E.coli	
Stem Cell Factor (KITLG)	S816-40N	Human		E.coli	
TFF1	T811-40N	Human		E.coli	
TFF3	T813-40N	Human		E.coli	
TGF Beta-3 (TGFB3)	T833-40N	Human		E.coli	
TGFB1	T831-40N	Human		E.coli	279-end
Thymosin-Beta 4 (TMSB4X)	T897-40N	Human		E.coli	
TNF	T860-40BN	Human		E.coli	
TNF-beta	T861-40BG	Mouse	GST	Sf9 insect cells using baculovirus	34-end
TNFRSF10B	T881-40BN	Human		E.coli	
TNFRSF1B	T863-40BN	Human		E.coli	
TSLP	T896-40N	Human		E.coli	
TWEAK (TNFSF12)	T893-40N	Human		E.coli	
TWEAK Receptor (TNFRSF12A)	T883-40N	Human		E.coli	
Uteroglobin (SCGB1A1)	U866-40N	Human		E.coli	
VEGF (121aa)	V810-40N	Human		E.coli	
VEGF (165aa)	V815-40N	Human		E.coli	
VEGFB	V811-40N	Human		E.coli	
WISP3	W810-40N	Human		E.coli	

# Affinity Proteins and Resins

Product Name	Cat# No
Ni-NTA Resin	NN01-RA091

**Compounds**



# Compounds

## Epigenetic Enzyme Activators

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
<b>BML-278</b>	S35-904C	N/F	C24H25NO4	>98% (TLC)	N-Benzyl-3,5-dicarbethoxy-4-phenyl-1,4-dihydropyridine
<b>CTPB</b>	P07-903	586976-24-1	C31H43ClF3NO2	98%	N-(4-chloro-3-trifluoromethyl-phenyl)-2-ethoxy-6-pentadecyl-benzamide
<b>Resveratrol</b>	S35-904	501-36-0	C14H12O3	>98%	3,4',5-Trihydroxy-trans-stilbene; 5-[(1E)-2-(4-Hydroxyphenyl)ethenyl]-1,3-benzenediol
<b>SirtAct</b>	S35-904E	839699-72-8	C20H25N5O2	98% (TLC)	2-Amino-N-cyclopentyl-1-(3-methoxypropyl)-1H-pyrrolo[2,3-b]quinoxaline-3-carboxamide; CAY10591

## Kinase Inhibitors

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
<b>6BIO</b>	G08-900	667463-62-9	C16H10BrN3O2	98% (TLC; 5% CH3OH3 in CH2Cl2, Rf=0.43); NMR (Conforms)	(2'Z,3'E)-6-Bromindirubin-3'-oxime
<b>7-CL-O-Nec-1</b>	R07-901B	852391-15-2	C13H12ClN3O	>98% (TLC); NMR (Conforms)	Nec-1s; 5-((7-Chloro-1H-indol-3-yl)methyl)-3-methylimidazolidine-2,4-dione
<b>A83-01</b>	A09-900	909910-43-6	C25H19N5S	>98% (TLC); NMR (Conforms)	3-(6-Methyl-2-pyridinyl)-N-phenyl-4-(4-quinolinyl)-1H-pyrazole-1-carbothioamide
<b>AAL-993</b>	F11-900C	269390-77-4	C20H16F3N3O	>98% (TLC); NMR (Conforms)	2-((4-Pyridyl)methyl)amino-N-(3 (trifluoromethyl)phenyl)benzamide
<b>AG-1296</b>	P12-900	146535-11-7	C16H14N2O2	>98% (TLC); NMR (Conforms)	6,7-Dimethoxy-3-phenylquinoxaline
<b>AG-1478</b>	E10-900B	175178-82-2	C16H14ClN3O2	>98% (TLC); NMR (Conforms)	"1. N-(3-Chlorophenyl)-6,7-dimethoxy-4-quinazolinanine 2. Free Base"
<b>AG-1517</b>	K11-900B	153436-54-5	C16H14BrN3O2 · HCl	>98% (TLC); NMR (Conforms)	"1. 4-(3-Bromoanilino)-6,7-dimethoxy-quinazoline hydrochloride 2. PD-153,035 3. SU-5271 4. ZM-252868"
<b>AG-490</b>	J02-901	133550-30-8	C17H14N2O3	>98% (TLC); NMR (Conforms)	(E)-2-Cyano-3-(3,4-dihydrophenyl)-N-(phenylmethyl)-2-propenamide
<b>AG-879</b>	N28-909B	148741-30-4	C18H24N2OS	>99% (TLC); NMR (Conforms)	alpha-Cyano-(3,5-di-t-butyl-4-hydroxy)thiocinnamide
<b>Alsterpaullone</b>	G09-901C	237430-03-4	C16H11N3O3	>96% (TLC); NMR (Conforms)	9-Nitro-7,12-dihydroindolo-[3,2-d][1]benzazepin-6(5)-one
<b>Amlexanox</b>	T02-901	68302-57-8	C16H14N2O4	>98% (TLC); NMR (Conforms)	2-Amino-7-(1-methylethyl)-5-oxo-5H-[1]benzopyrano[2,3-b]pyridine-3-carboxylic acid
<b>Apigenin</b>	K10-900C	520-36-5	C15H10O5	>98% (TLC); NMR (Conforms)	4',5,7-Trihydroxyflavone
<b>AR-A014418</b>	G09-901	487021-52-3	C12H12N4O4S	>98% (TLC); NMR (Conforms)	N-[(4-Methoxyphenyl) methyl]-N'-(5-nitro-2-thiazolyl) urea

# Kinase Inhibitors

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
AS-041164	P29-901	6318-41-8	C11H7NO4S	>98% (TLC); NMR (Conforms)	"1. 5-Benzo[1,3]dioxol-5-ylmethylene-thiazolidine-2,4-dione 2. NSC31098"
AS-605240	P27-901	648450-29-7	C12H7N3O2S	>98% (TLC); NMR (conforms)	5-(6-Quinoxalinylmethylene)-2,4-thiazolidine-2,4-dione
BAY 11-7082	C51-900	19542-67-7	C10H9NO2S	>98% (TLC); NMR (Conforms)	(E)-3-[(4-methylphenylsulfonyl)-2-propenenitrile
BML-259	C33-901	267654-00-2	C14H16N2OS	>98% (TLC: 5% CH3OH/CH2Cl2; Rf=0.46); NMR (Conforms)	N-(5-Isopropylthiazol-2-yl) phenylacetamide
BML-277	C48-901	516480-79-8	C20H14ClN3O2	≥98% (TLC; 1:1 ethylacetate: hexane; Rf=0.17); NMR (Conforms)	"1. 2-(4-(4-chlorophenoxy)phenyl)-1H-benzimidazole-5-carboxamide 2. Chk2 Inhibitor II"
BRD-7389	R15-901	376382-11-5	C24H18N2O2	>98% (TLC); NMR (Conforms)	1-[(2-Phenylethyl)amino]-3H-naphtho[1,2,3-de]quinoline-2,7-dione
CHIR-99021	G09-901B	252917-06-9	C22H18Cl2N8	>98% (TLC); NMR (Conforms)	6-[[2-[[4-(2,4-Dichlorophenyl)-5-(5-methyl-1H-imidazol-2-yl)-2-pyrimidinyl] amino] ethyl] amino]-3-pyridinecarbonitrile
CID-2011756	P72-901	638156-11-3	C22H21ClN2O3	≥98% (TLC); NMR (Conforms)	5-(3-Chlorophenyl)-N-[4-(4-morpholinylmethyl) phenyl]-2-furancarboxamide
CZC-54252 HCl	L10-901	1199911-27-9	C22H25ClN6O4S · HCl	≥98% (TLC); NMR (Conforms)	N-[-[[5-Chloro-2-[[2-methoxy-4-(4-morpholinyl) phenyl] amino]-4-pyrimidinyl] amino] phenyl] methanesulfonamide hydrochloride
Dabrafenib	B08-901	1195765-45-7	C23H20F3N5O2S2	>98% (TLC); NMR (conforms)	N-[3-[5-(2-amino-4-pyrimidinyl)-2-(1,1-dimethylethyl)-4-thiazolyl]-2-fluorophenyl]-2,6-difluoro-benzenesulfonamide; GSK2118436
Dasatinib	S19-901B	302962-49-8	C22H26ClN7O2S	>98% (TLC); NMR (Conforms)	"1. N-(2-chloro-6-methylphneyl)-2-(6-(4-(2-hydroxyethyl)-piperazin-1-yl)-2-methylpyrimidin-4-ylamino) thiazole-5-carboxamide; 2. BMS-354825"
Deguelin	A16-900B	522-17-8	C23H22O6	>98% (TLC); NMR (Conforms)	None
Dibutyryl cAMP	P51-901	16980-89-5	C18H23N5O8P · Na	>98% (HPLC); NMR (conforms)	N6, O2'-Dibutyryl adenosine 3', 5'-cyclic monophosphate sodium salt; Bucladesine Na
Dimethylsphingosine	S17-900	119567-63-4	C20H41NO2	>98% (TLC); NMR (Conforms)	(2S,3R,E)-2-(dimethylamino)octadeca-4-ene-1,3,diol
DMPQ 2 HCl	P13-901	137206-97-4	C16H14N2O2 · 2HCl	>98% (TLC); NMR (Conforms)	5,7-Dimethoxy-3-(4-pyridinyl) quinoline dihydrochloride
Doramapimod	M39-900	285983-48-4	C31H37N5O3	>98% (TLC); NMR (Conforms)	"1. 1-[5-tert-butyl-2-(4-methylphenyl) pyrazol-3-yl]-3-[4-(2-morpholin-4-ylethoxy) naphthalen-1-yl] urea 2. BIRB-796"
Dorsomorphin	P47-900	866405-64-3	C24H25N5O	>98% (TLC); NMR (Conforms)	"1. 6-[4-[2-(1-Piperidinyl) ethoxy] phenyl]-3-(4-pyridinyl)-pyrazolo[1,5-a] pyrimidine 2. BML-275"
Erlotinib HCl	E10-900C	183319-69-9	C22H23N3O4 · HCl	>98% (TLC); NMR (Conforms)	N-(3-ethynylphenyl)-6,7-bis(2-methoxyethoxy)quinazolin-4-amine, hydrochloride
Everolimus	F17-901D	159351-69-6	C53H83NO14	≥98% (HPLC); NMR (Conforms)	40-O-(2-hydroxyethyl)rapamycin)

# Kinase Inhibitors

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
Fasudil HCl	R10-900	105628-07-7	C14H17N3O2S · HCl	>98% (TLC); NMR (Conforms)	"1. 1-(5-Isoquinolinesulfonyl) homopiperazine HCl 2. HA-1077"
Flavopiridol HCl	C22-900	131740-09-5	C21H20ClNO5 · HCl	>98% (HPLC); NMR (Conforms)	2-(2-Chlorophenyl)-5,7-dihydroxy-8-[(3S, 4R)-3-hydroxy-1-methyl-4-piperidinyl]-4H-1-benzopyran-4-one, hydrochloride; L 86-8275; Alvocidib; NSC-649890
Gefitinib	E10-900	184475-35-2	C22H24ClFN4O3	>98% (TLC); NMR (Conforms)	"1. N-(3-chloro-4-fluorophenyl)-7-methoxy-6-(3-morpholinopropoxy) quinazolin-4-amine; 2. ZD1839"
Genistein	K11-900	446-72-0	C15H10O5	98% (TLC); NMR (Conforms)	"1. 5,7-Dihydroxy-3-(4-hydroxyphenyl)-4H-1-benzopyran-4-one 2. 4',5,7-Trihydroxyisoflavone"
GF-109203X	P61-900	133052-90-1	C25H24N4O2	98% (TLC); NMR (Conforms)	"1. 2-[1-(3-Dimethylaminopropyl)-1H-indol-3-yl]-3-(1H-indol-3-yl) maleimide 2. Bisindolylmaleimide I"
GSK-690693 HCl	A16-900	937174-76-0	C21H27N7O3 · HCl	99% (HPLC); NMR (Conforms)	(S)-4-(2-(4-amino-1,2,5-oxadiazol-3-yl)-1-ethyl-7-(piperidin-3-ylmethoxy)-1H-imidazo[4,5-c]pyridin-4-yl)-2-methylbut-3-yn-2-ol. HCl
GW-441756	N28-909	504433-23-2	C17H13N3O	98% (TLC; 5% CH3OH: CH2Cl2; Rf=0.4) mixture of E and Z isomers	1,3-Dihydro-3-[(1-methyl-1H-indol-3-yl)methylene]-2H-pyrrolo[3,2-b]pyridin-2-one
GW-5074	R01-901B	220904-83-6	C15H8Br2INO2	98% (TLC); NMR (Conforms)	3-(3,5-Dibromo-4-hydroxy-benzylidene)-5-iodo-1,3-dihydro-indol-2-one
GW2580	C74-901	870483-87-7	C20H22N4O3	≥98% (TLC); NMR (Conforms)	5-(3-methoxy-4-((4-methoxybenzyl)oxy)benzyl)pyrimidine-2,4-diamine
H-89 2HCl	P51-900B	127243-85-0	C20H20BrN3O2S · 2HCl	>98% (TLC); NMR (Conforms)	N-[2-[[3-(4-Bromophenyl)-2-propenyl]amino]ethyl]-5-isoquinolinesulfonamide dihydrochloride
HA-100 2HCl	K10-900F	84468-24-6	C13H15N3O2S · 2HCl	>98% (HPLC); NMR (conforms)	5-(1-Piperazinylsulfonyl)isoquinoline dihydrochloride
Hypericin	P61-900C	548-04-9	C30H16O8	>98% (TLC); NMR (Conforms)	1,3,4,6,8,13-Hexahydroxy-10,11-dimethylphenanthro [1,10,9,8-opqra] perylene-7,14-dione
IM-12	G09-901D	1129669-05-1	C22H20FN3O2	>98% (HPLC); NMR (conforms)	3-(4-Fluorophenylethylamino)-1-methyl-4-(2-methyl-1H-indol-3-yl)-1H-pyrrole-2,5-dione
Imatinib mesylate	A03-901B	220127-57-1	C29H31N7O · CH3SO3H	98% (TLC); NMR (Conforms)	"1. 4-[(4-Methyl-1-piperazinyl) methyl]-N-[4-methyl-3-[[4-(3-pyridinyl)-2-pyrimidinyl] amino]-phenyl] benzamide methane-sulfonate 2. STI-571"
Indirubin-3'-monoxime	C29-901	160807-49-8	C16H11N3O2	>98% (TLC); NMR (Conforms)	Indirubin-3'-oxime 3-[1,3-Dihydro-3-(hydroxyimino)-2H-indol-2-ylidene]-1,3-dihydro-2H-indol-2-one
IQ-1S	M35-901B	1421610-21-0	269.2	>98% (TLC); NMR (Conforms)	11H-Indeno[1,2-b]quinoxalin-11-one oxime, sodium salt
JNK Inhibitor IX	M34-901	312917-14-9	C20H16N2OS	>98% (TLC); NMR (Conforms)	N-(3-cyano-4,5,6,7-tetrahydrobenzo[b]thien-2-yl)-1-naphthalenecarboxamide
JX-401	M39-901	349087-34-9	C21H25NO2S	>98% (TLC); NMR (Conforms)	1-[2-Methoxy-4-(methylthio) benzoyl]-4-(phenylmethyl) piperidine
K252A	K10-900D	99533-80-9	C27H21N3O5	>98% (TLC); NMR (Conforms)	None

# Kinase Inhibitors

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
<b>K252B</b>	K10-900E	99570-78-2	C26H19N3O5	>98% (TLC); NMR (Conforms)	None
<b>KT-5823</b>	P78-900	126643-37-6	C29H25N3O5	>98% (TLC); NMR (Conforms)	None
<b>KT5720</b>	P51-900	108068-98-0	C32H31N3O5	>98% (TLC); NMR (Conforms)	None
<b>KU-55933</b>	A26-901	587871-26-9	C21H17N3O3S2	>98% (TLC); NMR (conforms)	2-(4-Morpholinyl)-6-(1-thianthrenyl)-4H-pyran-4-one
<b>Lapatinib</b>	E10-900D	231277-92-2	C29H26ClFN4O4S	98% (TLC); NMR (Conforms)	"1. N-[3-chloro-4-[(3-fluorophenyl) methoxy]phenyl]-6-[5-[(2-methylsulfonyl-ethylamino)methyl]-2-furyl]quinazolin-4-amine 2. GW-572016"
<b>Lenvatinib</b>	K01-901	417716-92-8	C21H19ClN4O4	>98% (TLC); NMR (conforms)	4-[3-Chloro-4-[[[(cyclopropylamino) carbonyl]amino]phenoxy]-7-methoxy-6-quinolinecarboxamide; E7080
<b>Lestaurtinib</b>	F12-901	111358-88-4	C26H21N3O4	>98% (TLC); NMR (Conforms)	"1. CEP-701 2. KT-5555"
<b>LFM-A13</b>	B10-901	62004-35-7	C11H8Br2N2O2	98% (TLC: 2:8 CH3OH:CH2Cl2; Rf=0.53); NMR (Conforms)	2-Cyano-N-(2,5-dibromophenyl)-3-hydroxy-2-butenamide
<b>LH-846</b>	C67-901	639052-78-1	C16H13ClN2OS	>98% (TLC); NMR (Conforms)	N-(5-Chloro-6-methyl-2-enzothiazolyl) benzeneacetamide
<b>Locostatin</b>	R01-901C	133812-16-5	C14H15NO3	>98% (TLC); NMR (conforms)	(4S)-3-[(E)-But-2-enoyl]-4-benzyl-2-oxazolidinone; UIC-1005
<b>LY-294002</b>	P27-900C	154447-36-6	C19H17NO3	>98% (TLC); NMR (Conforms)	2-(4-Morpholinyl)-8-phenyl-4H-1-benzopyran-4-one
<b>LY-364947</b>	T07-901	396129-53-6	C17H12N4	>98% (TLC); NMR (Conforms)	4-(3-Pyridin-2-yl-1H-pyrazol-4-yl)quinoline; HTS-466284
<b>LYS6K2</b>	R15-900	NF	C21H19F4N7	>97% (HPLC); NMR (Conforms)	4-(4-(4-(4-fluoro-3-(trifluoromethyl) phenyl)-1-methyl-1H-imidazol-2-yl) piperidin-1-yl)-1H-pyrazolo [3,4-d] pyrimidine
<b>MLS-573151</b>	C08-909	NF	C21H19N3O2S	>98% (TLC); NMR (Conforms)	4-[3,5-diphenyl-4,5-dihydropyrazol-1-yl] benzenesulfonamide
<b>Necrostatin-1</b>	R07-901	4311-88-0	C13H13N3OS	>98% (TLC: 1:9, CH3OH:CH2Cl2; Rf=0.6); NMR (Conforms)	"1. 5-(1H-Indol-3-ylmethyl)-3-methyl-2-thioxo-4-imidazolidinone 2. Methylthiohydantoin-DL-tryptophan"
<b>Nilotinib</b>	A03-901	641571-10-0	C28H22F3N7O	>98% (TLC); NMR (Conforms)	"1. 4-methyl-N-[3-(4-methyl-1H-imidazol-1-yl)- 5-(trifluoromethyl) phenyl]-3-[(4-pyridin-3-ylpyrimidin-2-yl) amino] benzamide 2. AMN-107"
<b>NPC-15437</b>	P61-900B	141774-20-1	C25H50N4O2 · 2HCl	98% (TLC); NMR (Conforms)	(S)-2,6-Diamino-N-[(1-(1-oxotridecyl)-2-piperidinyl) methyl] hexanamide dihydrochloride hydrate
<b>NVP-BEZ235</b>	F17-900	915019-65-7	C30H23N5O	>98% (TLC); NMR (Conforms)	2-Methyl-2-[4-[3-methyl-2-oxo-8-(quinolin-3-yl)-2,3-dihydro-1H-imidazo[4,5-c] quinolin-1-yl] phenyl] propanenitrile
<b>Parthenolide</b>	103-901B	20554-84-1	C15H20O3	>98% (TLC); NMR (conforms)	Sesquiterpene lactone found in feverfew (Tanacetum parthenium)

# Kinase Inhibitors

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
<b>PD-325901</b>	M02-901B	391210-10-9	C16H14F3IN2O4	≥98% (HPLC); NMR (Conforms)	N-[(2R)-2,3-Dihydroxypropoxy]-3,4-difluoro-2-[(2-fluoro-4-iodophenyl)amino]-benzamide
<b>PD-98059</b>	M02-900	167869-21-8	C16H13NO3	>98% (TLC); NMR (Conforms)	2-(2-Amino-3-methoxyphenyl)-4H-1-benzopyran-4-one; 2'-Amino-3'-methoxyflavone
<b>Phorbol-12-myristate-13 acetate</b>	P61-900D	16561-29-8	C36H56O8	>99% (HPLC); NMR (conforms)	12-O-Tetradecanoylphorbol-13-acetate; PMA; TPA
<b>Piceatannol</b>	I01-901	10083-24-6	C14H12O4	>98% (TLC); NMR (Conforms)	3,3',4,5'-Tetrahydroxy-trans-stilbene
<b>Picropodophyllotoxin</b>	I02-901	477-47-4	C22H22O8	>97% (TLC); NMR (Conforms)	(5R,5aS,8aR,9R)-5,8,8a,9-Tetrahydro-9-hydroxy-5-(3,4,5-trimethoxyphenyl)-furo[3',4':6,7]naphtho[2,3-d]-1,3-dioxol-6(5aH)-one
<b>PKC-412</b>	K10-900B	120685-11-2	C35H30N4O4	>98% (TLC); NMR (Conforms)	"1. 4'-N-Benzoyl staurosporine 2. Midostaurin 3. CGP-41251"
<b>PP1</b>	S19-900	172889-26-8	C16H19N5	98% (TLC); NMR (Conforms)	4-Amino-1-tert-butyl-3-(4-methylphenyl)pyrazolo [3,4-d] pyrimidine
<b>PP2</b>	S19-900B	172889-27-9	C15H16ClN5	>98% (TLC); NMR (Conforms)	3-(4-chlorophenyl)1-(1,1-dimethylethyl)-1H-pyrazolo [3,4-d] pyrimidin-4-amine
<b>PX-866</b>	P27-900	502632-66-8	C29H35NO8	>98% (TLC); NMR (Conforms)	(1E,4S,4aR,5R,6aS,9aR)-5-(acetyloxy)-1-[(di-2-propen-1-ylamino)methylene]-4,4a,5,6,6a,8,9,9-a-octa-hydro-11-hydroxy-4-(methoxymethyl)-4a,6a-dimethyl-cyclopenta [5,6] naphtha [1,2-c] pyran-2,7,10(1H)-trione
<b>R-59022</b>	D21-900	93076-89-2	C27H26FN3OS	>98% (HPLC); NMR (conforms)	6-[2-[4-[(4-Fluorophenyl)phenylmethylene]-1-piperidinyl]ethyl]-7-methyl-5H-thiazolo[3,2-a]pyrimidin-5-one
<b>Rapamycin</b>	F17-901	53123-88-9	C51H79NO13	>98% (TLC: 10% CH3OH in CH2Cl2; Rf=0.51); NMR (Conforms)	Sirolimus; Fermentation product
<b>Regorafenib</b>	K01-900	755037-03-7	C21H15ClF4N4O3	>98% (HPLC,TLC); NMR (conforms)	4-[4-[[[4-chloro-3-(trifluoromethyl)phenyl]amino]carbonyl]amino]-3-fluorophenoxy]-N-methyl-2-pyridinecarboxamide; BAY 73-4506
<b>Reversine</b>	A28-900	656820-32-5	C21H27N7O	>98% (HPLC, TLC); NMR (conforms)	2-(4-Morpholinoanilino)-6-cyclohexylaminopurine
<b>Ro 31-6045</b>	R21-901	113963-68-1	C21H15N3O2	>98% (HPLC, TLC); NMR (conforms)	1-Methyl-3,4-bis(indol-3-yl)maleimide; Bisindolylmaleimide V; BIM V
<b>Roscovitine</b>	C22-901	186692-46-6	C19H26N6O	>99% (TLC); NMR (Conforms)	(R)-2-((6-(benzylamino)-9-isopropyl-9H-purin-2-yl) amino) butan-1-ol
<b>Ruxolitinib</b>	J01-901	941678-49-5	C17H18N6	>98% (HPLC,TLC); NMR (conforms)	(3R)-3-cyclopentyl-3-[4-(7H-pyrrolo[2,3-d]pyrimidin-4-yl)pyrazol-1-yl]propanenitrile
<b>SB-202190</b>	M39-900B	152121-30-7	C20H14N3OF	>98% (TLC: 10% CH3OH in CH2Cl2, Rf=0.25); NMR (Conforms)	4-[4-(4-Fluorophenyl)-5-(4-pyridinyl)-1H-imidazol-2-yl] phenol
<b>SB-203580</b>	M39-900C	152121-47-6	C21H16N3OSF	>98% (TLC); NMR (Conforms)	4-(4-fluorophenyl)-2-(4-methylsulfinylphenyl)-5-(4-pyridyl)imidazole

# Kinase Inhibitors

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
SC-514	I03-901	354812-17-2	C9H8N2OS2	>98% (TLC; 5% CH3OH in CH2Cl2; Rf=0.27); NMR (Conforms)	5-(Thien-3-yl)-3-aminothiophene-2-carboxamide
Sorafenib	R01-900	284461-73-0	C21H16ClF3N4O3	>99% (TLC); NMR (Conforms)	"1. 4-(4-(3-(4-chloro-3 (trifluoromethyl) phenyl) ureido) phenoxy)-N-methylpicolinamide 2. BAY 43-9006"
SP-600125	M33-900	129-56-6	C14H8N2O	>98% (TLC); NMR (Conforms)	"1. Anthra [1-9-cd] pyrazol-6(2H)-one 2. 1,9-Pyrazoloanthrone"
SR-3576	M35-901	1164153-22-3	C27H27N5O5	>98% (TLC); NMR (Conforms)	3-(5-(3-m-Tolylureido)-1H-indazol-1-yl)-N-(3,4,5-trimethoxyphenyl)benzamide
Src I1	S19-901	179248-59-0	C22H19N3O3	≥98% (TLC; 5% CH3OH in CH2Cl2; Rf=0.43); NMR (Conforms)	6,7-Dimethoxy-N-(4-phenoxyphenyl)-4-quinazolinamine
SRPIN340	S21-901	218156-96-8	C18H18F3N3O	>98% (TLC); NMR (Conforms)	N-[2-(1-Piperidiny)-5-(trifluoromethyl) phenyl]-4-pyridinecarboxamide
Staurosporine	K10-900	62996-74-1	C28H26N4O3	>98% (TLC); NMR (Conforms)	
STO-609	C17-900	52029-86-4	C19H10N2O3 · CH3COOH	98% (TLC); NMR (Conforms)	7-oxo-7H-benzimidazo[2,1-a]benz[de]isoquinoline-3-carboxylic acid acetate
SU-1498	F11-900E	NF	C25H30N2O2	>98% (TLC); NMR (Conforms)	(E)-N-(3-Phenylpropyl)-alpha-cyano-3',5'-diisopropyl-4'-hydroxycinnamamide
SU-4312	F11-900F	5812-07-07	C17H16N2O	>98% (TLC); NMR (Conforms)	"1. 3-(4-Dimethylaminobenzylidene)-1,3-dihydroindol-2-one 2. DMBI"
Sunitinib malate	F11-900G	341031-54-7	C22H27FN4O2 · C4H6O5	>98% (TLC); NMR (Conforms)	"1. N-[2-(Diethylamino)ethyl]-5-[(Z)-(5-fluoro-1,2-dihydro-2-oxo-3H-indol-3-ylidene) methyl]-2,4-dimethyl-1H-pyrrole-3-carboxamide malate 2. SU-11248"
TG003	C57-901	719277-26-6	C13H15NO2S	>98% (HPLC); NMR (Conforms)	(Z)-1-(3-Ethyl-5-methoxy-2,3-dihydrobenzothiazol-2-ylidene)propan-2-one
U-0126	M02-901	109511-58-2	C18H16N6S2	>98% (TLC); NMR (Conforms)	1,4-Diamino-2,3-dicyano-1,4-bis [2-aminophenylthio] butadiene
Vandetanib	F11-900	443913-73-3	C22H24BrFN4O2	>98% (TLC); NMR (Conforms)	"1. 4-(4-Bromo-2-fluoroanilino)-6-methoxy-7-[(1-methylpiperidin-4-yl) methoxy] quinazoline 2. ZD6474"
Vatalanib 2HCl	F11-900D	212141-51-0	C20H15ClN4·2HCl	>98% (TLC); NMR (Conforms)	"1. N-(4-Chlorophenyl)-4-(4-pyridinylmethyl)-1-phthalazinamine dihydrochloride 2. CGP-79787 3. PTK-787 4. ZK222584"
WHI-P131	J03-901	202475-60-3	C16H15N3O3 · HCl	>98% (TLC); NMR (Conforms)	"1. 4-[(4'-hydroxyphenyl)amino]-6,7-dimethoxyquinazoline, Hydrochloride 2. Janex-1"
Wortmannin	P27-900B	19545-26-7	C23H24O8	>98% (TLC); NMR (Conforms)	None
Y-27632 2HCl	R10-900B	129830-38-2	C14H21N3O · 2HCl	99% (HPLC); NMR (Conforms)	(R)-(+)-trans-4-(1-Aminoethyl)-N-(-4-pyridyl) cyclohexanecarboxamide dihydrochloride

# Kinase Inhibitors

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
<b>ZM-306416</b>	F11-900B	690206-97-4	C16H13ClFN3O2	>98% (TLC); NMR (Conforms)	4-[(4'-Chloro-2'-fluoro) phenylamino]-6,7-dimethoxyquinazoline
<b>ZM-336372</b>	R01-901	331771-20-1	C23H23N3O3	98% (TLC); NMR (Conforms)	3-(Dimethylamino)-N-[3-[(4-hydroxybenzoyl)-amino]-4-methylphenyl]-benzamide
<b>ZM-447439</b>	A31-901	331771-20-1	C29H31N5O4	>98% (TLC); NMR (Conforms)	(N-[4-[[6-Methoxy-7-[3-(4-morpholinyl)propoxy]-4-quinazolinyl] amino] phenyl] benzamide

# Phosphatase Inhibitors

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
<b>Cantharidin</b>	P16-902C	56-25-7	C10H12O4	>98% (HPLC)	3a,7a-Dimethylhexahydro-4,7-epoxyisobenzofuran-1,3-dione
<b>NSC-663284</b>	C04-902	383907-43-5	C15H16ClN3O3	>98%	6-Chloro-7-[[2-(4-morpholinyl)ethyl] amino]-5,8-quinolinedione
<b>Okadaic acid, free acid</b>	P16-902	78111-17-8	C44H68O13	>98% (TLC)	None

# Epigenetic Enzyme Inhibitors

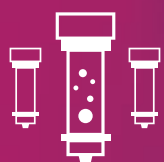
Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
<b>2'-Deoxy-5-fluoro-cytidine</b>	D353-904D	10356-76-0	C9H12FN3O4	>98%	5-Fluoro-2'-deoxycytidine; FCDR; FdCyd; NSC-48006
<b>5-Azacytidine</b>	D353-904	320-67-2	C8H12N4O5	>98% (TLC)	4-Amino-1-(beta-D-ribofuranosyl)-1,3,5-triazin-2(1H)-one, Ladakamycin
<b>AK-7</b>	S36-904	420831-40-9	C19H21BrN2O3S	98% (TLC)	N-(3-Bromophenyl)-3-[(hexahydro-1H-azepin-1-yl)sulfonyl]benzamide
<b>APICIDIN</b>	H83-904B	183506-66-3	C34H49N5O6	≥98%	Cyclo[(2S)-2-Amino-8-oxodecanoyl-1-methoxy-L-tryptophyl-L-isoleucyl-(2R)-2-piperidinecarbonyl]
<b>BIX-01294 3HCl</b>	S344-903	935693-62-2	C28H38N6O2.3HCl	>98%	2-(Hexahydro-4-methyl-1H-1,4-diazepin-1-yl)-6,7-dimethoxy-N-[1-(phenylmethyl)-4-piperidinyl]-4-quinazolinamine trihydrochloride
<b>BML-210</b>	H83-904C	537034-17-6	C20H25N3O2	98%	N1-(2-aminophenyl)-N8-phenyloctanedi-amide
<b>C-646</b>	K311-903	328968-36-1	C24H19N3O6	>98%	4-[4-[[5-(4,5-Dimethyl-2-nitrophenyl)-2-furanyl]methylene]-4,5-dihydro-3-methyl-5-oxo-1H-pyrazol-1-yl]benzoic acid
<b>Chaetocin</b>	K421-904C	28097-03-2	C30H28N6O6S4	>99%	Lys methyltransferase inhibitor
<b>CPH2</b>	K311-903B	357649-93-5	C14H14ClN3S	>98% (TLC)	Cyclopentylidene-[4-(4'-chlorophenyl)thiazol-2-yl]hydrazone
<b>CUDC-101</b>	H83-904	1012054-59-9	C24H26N4O4	>99% (HPLC)	7-((4-((3-ethynylphenyl)amino)-7-methoxyquinazolin-6-yl)oxy)-N-hydroxyheptanamide
<b>Daminozide</b>	K421-904B	1596-84-5	C6H12N2O3	98%	1-(2,2-Dimethylhydrazide)butanedioic acid
<b>Decitabine</b>	D353-904B	2353-33-5	C8H12N4O4	>98%	5-Aza-2'-deoxycytidine; NSC-127716
<b>DMOG</b>	J48-904	89464-63-1	C6H9NO5	>98%	N-(2-Methoxy-2-oxoacetyl)glycine methyl ester; Dimethyloxalylglycine

# Epigenetic Enzyme Inhibitors

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
EX-527	S35-904F	49843-98-3	C13H13ClN2O	98% (TLC)	6-Chloro-2,3,4,9-tetrahydro-1H-carbazole-1-carboxamide (racemic)
Garcinol	K311-903C	78824-30-3	C38H50O6	>97% (TLC)	Camboginol, 3-(3,4-dihydroxybenzoyl)-4-hydroxy-8,8-dimethyl-1,7-bis(3-methyl-2-buten-1-yl)-5-[(2S)-5-methyl-2-(1-methylethenyl)-4-hexen-1-yl]-bicyclo[3.3.1]non-3-ene-2,9-dione
GSK-J1 (Free acid)	J48-904B	1373422-53-7	C22H23N5O2	99%	N-[2-(2-Pyridinyl)-6-(1,2,4,5-tetrahydro-3H-3-benzazepin-3-yl)-4-pyrimidinyl]-β-alanine
GSK-J4 (ethylester)	J48-904C	1373423-53-0	C24H27N5O2	>99%	N-[2-(2-Pyridinyl)-6-(1,2,4,5-tetrahydro-3H-3-benzazepin-3-yl)-4-pyrimidinyl]-β-alanine ethyl ester
HPA (HEXYL-4-PENTYNOIC ACID)	H83-904D	96017-59-3	C11H18O2	>97% (TLC)	2-(2-propynyl)octanoic acid (racemic); HPA
IOX1 (8H5QCA)	J48-904D	5852-78-8	C10H7NO3	>98% (HPLC)	5-Carboxy-8-hydroxyquinoline, 8-Hydroxy-5-quinolinecarboxylic acid
ML-324	J48-904E	1222800-79-4	C21H23N3O2	>98% (TLC)	N-(3-(dimethylamino)propyl)-4-(8-hydroxyquinolin-6-yl)benzamide
MS-275 (Entinostat)	H83-904E	209783-80-2	C21H20N4O3	>98% (TLC)	Pyridin-3-ylmethyl 4-((2-aminophenyl)carbamoyl)benzylcarbamate; Entinostat
Panobinostat	H83-904G	404950-80-7	C21H23N3O2	>98% (TLC)	(E)-N-hydroxy-3-(4-((2-(2-methyl-1H-indol-3-yl)ethylamino)methyl)phenyl) crylamide; LBH-589
Phenylbutyrate NA	H83-904H	1716-12-7	C10H11NaO2	>98% (TLC; 5% CH3OH in CH2Cl2, Rf=0.36)	4-Phenylbutyric acid, sodium salt
RG-108	D353-904C	48208-26-0	C19H14N2O4	>98% (TLC)	N-Phthalyl-L-tryptophan
Romidepsin (FK228)	H83-904I	128517-07-7	C24H36N4O6S2	>98% (HPLC)	Cyclo[(2Z)-2-amino-2-butenoyl-L-valyl-(3S,4E)-3-hydroxy-7-mercapto-4-heptenoyl-D-valyl-D-cysteinyl], cyclic (3-5) disulfide
Sirtinol	S35-904B	410536-97-9	C26H22N2O2	>98% (TLC)	2-[[[(2-Hydroxy-1-naphthalenyl)methylene]amino]-N-(1-phenylethyl)benzamide
Splitomicin	S35-904G	5690-03-09	C13H10O2	>98% (TLC)	1,2-Dihydro-3H-naphtho[2,1-b]pyran-3-one
TC-E 5003	P3695-903	17328-16-4	C16H14Cl2N2O4S	>98% (TLC)	N,N'-(Sulfonyldi-4,1-phenylene)bis(2-chloroacetamide)
Tranylcypromine	K421-904	13492-01-8	C9H11N · 1/2H2SO4	>98% (TLC)	Trans (±)-2-phenylcyclopropanamine; 2-PCPA
Trichostatin A	H83-904J	58880-19-6	C17H22N2O3	>99%	(2E,4E,6R)-7-(4-(Dimethylamino)phenyl)-N-hydroxy-4,6-dimethyl-7-oxo-2,4-heptadienamide
Tubastatin A · HCL	H83-904K	1239262-52-2	C20H21N3O2·HCl	99% (HPLC)	N-hydroxy-4-((2-methyl-1,2,3,4-tetrahydropyrido[4,3-b]indol-5-yl)methyl) benzamide hydrochloride
UNC1999	D353-904E	1431612-23-5	C33H43N7O2	>98% (HPLC)	N-[(6-methyl-2-oxo-4-propyl-1H-pyridin-3-yl)methyl]-1-propan-2-yl-6-[6-(4-propan-2-ylpiperazin-1-yl)pyridin-3-yl]indazole-4-carboxamide
Valproate NA	H83-904L	1069-66-5	C8H15NaO2	>98% (TLC)	Sodium 2-propylpentanoate
Vorinostat (SAHA)	H83-904M	149647-78-9	C14H20N2O3	>98%	N-Hydroxy-N'-phenyloctanediamide

# Kinase Activators

Product Name	Cat# No	CAS Registration #	Molecular Formula	Purity	Chemical Name
<b>8-pCPT-2-O-Me-cAMP-AM</b>	R46-905	1152197-23-3	C20H21ClN5O8PS	>97% (HPLC)	8-(4-Chlorophenylthio)-2'-O-methyl-adenosine-3', 5'-cyclic monophosphate acetoxymethyl ester
<b>AICAR</b>	P47-900B	2627-69-2	C9H14N4O5	>99% (TLC); NMR (Conforms)	"1. 5-Aminoimidazole-4-carboxamide 1-β-D-ribofuranoside 2. Acadesine"
<b>MHY-1485</b>	F17-901B	326914-06-1	C17H21N7O4	>98% (TLC); NMR (conforms)	4,6-Di-4-morpholinyl-N-(4-nitrophenyl)-1,3,5-triazin-2-amine
<b>SC-79</b>	A16-901	305834-79-1	C17H17ClN2O5	98% (TLC); NMR (Conforms)	2-Amino-6-chloro-α-cyano-3-(ethoxycarbonyl)-4H-1-benzopyran-4-acetic acid, ethyl ester



# Reagents

# Reagents

## Assay Buffer and Cofactors

Product Name	Cat# No	Application
Acetyltransferase Assay Buffer	A01-09	Acetyltransferase Assay
Acetyltransferase Dilution Buffer	A21-09	Acetyltransferase Assay
AMP Solution	A46-09	Kinase Assay
ATP Stock Solution	A50-09	Kinase Assay
Ca <sup>2+</sup> /Calmodulin Solution II, 10x	C02-39B	Kinase Assay
CaCl <sub>2</sub> Solution	C02-09	Enzyme Assay
cAMP Solution	A47-09	Enzyme Assay
cGMP Solution	G47-09	Enzyme Assay
DTT Solution	D86-09B	
DTT Solution	D86-09	
DUB Reaction Buffer (4X)	D01-09	
Fructose Kinase Assay Buffer I	F01-09	Enzyme Assay
Fructose Kinase Dilution Buffer I	F21-09	Enzyme Assay
GTP Solution	G50-09	Enzyme Assay
GTP Solution II	G50-09B	Enzyme Assay
Kinase Assay Buffer I	K01-09	Kinase Assay
Kinase Assay Buffer II	K02-09	Kinase Assay
Kinase Assay Buffer III	K03-09	
Kinase Dilution Buffer I	K21-09	Kinase Assay
Kinase Dilution Buffer II	K22-09	Kinase Assay
Name	Catalog Number	Application
Kinase Dilution Buffer III	K23-09	Kinase Assay
Kinase Dilution Buffer IV	K24-09	Kinase Assay
Kinase Dilution Buffer IX	K29-09	
Kinase Dilution Buffer V	K25-09	Kinase Assay
Kinase Dilution Buffer VI	K26-09	Kinase Assay
Kinase Dilution Buffer VII	K27-09	Kinase Assay
Kinase Dilution Buffer VIII	K28-09	Kinase Assay
Kinase Dilution Buffer X	K20-09	Kinase Assay
Lipid Dilution Buffer	L21-09	
Lipid Dilution Buffer II	L22-09	
Lipid Kinase Buffer	L01-09	
Methyltransferase Assay Buffer I	M01-09	Kinase Assay
Methyltransferase Assay Buffer II	M02-09	Kinase Assay
Methyltransferase Assay Buffer III	M03-09	Kinase Assay
Methyltransferase Assay Buffer IV	M04-09	
Methyltransferase Dilution Buffer I	M21-09	Kinase Assay
Methyltransferase Dilution Buffer II	M22-09	Kinase Assay
Methyltransferase Dilution Buffer III	M23-09	Kinase Assay
Methyltransferase Dilution Buffer IV	M24-09	
METTL12 Substrate	M332-58	
MnCl <sub>2</sub> Solution	M40-09	Enzyme Assay

# Assay Buffer and Cofactors

Product Name	Cat# No	Application
Neddylation Buffer	N01-09	
Phosphatase Assay Buffer I	P01-09	Phosphatase Assay
Phosphatase Assay Buffer II	P02-09	Phosphatase Assay
Phosphatase Assay Buffer III	P03-09	Phosphatase Assay
Phosphatase Dilution Buffer I	P21-09	Phosphatase Assay
Phosphatase Dilution Buffer II	P22-09	Phosphatase Assay
Phosphatase Dilution Buffer III	P23-09	Phosphatase Assay
PKC Lipid Activator, 10x	L51-39	Kinase Assay

# Assay Substrates

## Lipid Substrates

Product Name	Cat# No	Application
DLG (Dilauroyl-sn-glycerol)	D430-59	
PI (Phosphatidylinositol)	P420-59	
PI(3)P	P421-59	
PI(3)P:PS	P426-59	
PI(4)P	P422-59	
PI(4)P:PS	P427-59	
PI(4,5)P2	P424-59	
PI(4,5)P2:PS	P429-59	
PI(5)P	P423-59	
PI(5)P:PS	P428-59	
PI:PS	P425-59B	
PS (Phosphatidylserine) Solution	L11-39	

## Oligo Substrates

Product Name	Cat# No	Application
DNMT Substrate-1	D53-58	DNA methyltransferase assay (50pmole per reaction).
METTL3/METTL14 Substrate	M323-58	A methyltransferase METTL3/METTL14 complex assay.
Poly (1:1 dl, dC) Acid	P62-58	DNA methyltransferase assay (0.25ug per reaction).

## Carbohydrate Substrates

Product Name	Cat# No	Application
D-Fructose-6-phosphate	F520-59	PFKFB Assay
D-Glucose-6-phosphate	F521-59	PFKFB Assay

# Peptide Substrates

Product Name	Cat# No	Application
3CLpro FRET Substrate	CL01-58	COVID-19 Assays
Abltide	A02-58	Kinase Assay
Acetylated Histone H3 Peptide I	H12-358	Acetylation/Methylation Assays
Acetylated Histone H4 Peptide I	H13-358	Acetylation/Methylation Assays
AKT (PKB) Substrate	A05-58	Kinase Assay
AKT (SGK) Substrate	A08-58	Kinase Assay
AMARA Peptide	A11-58	Kinase Assay
Amyloid Beta Peptide (1-42)	A42-58	
Autocamtide 2	A15-58	Kinase Assay
Axltide	A16-58	Kinase Assay
CATCHtide	C02-58	Kinase Assay
CDKtide	C06-58	Kinase Assay
Chktide	C10-58	Kinase Assay
CK1tide	C07-58	Assay
CK2 Substrate	C08-58	Kinase Assay
CREBtide	C50-58	Kinase Assay
Crosstide	C51-58	Kinase Assay
CSKtide	C63-58	Kinase Assay
DYRKtide	D96-58	Kinase Assay
EF2tide	E01-58	Kinase Assay
EIF2S Peptide	E23-58	Kinase Assay
GRKtide	G46-58	Kinase Assay
GS Peptide	G60-58	Kinase Assay
GSK3 Substrate	G50-58	Kinase Assay
HER2 Substrate	E27-58	Kinase Assay
Histone H1 Peptide	H10-58	Assay
Histone H3 Peptide (1-21)	H12-58	Acetylation/Methylation Assays
Histone H4 Peptide (1-21)	H13-58	Acetylation/Methylation Assays
HSP27tide	H31-58	Kinase Assay
IGF1Rtide	I15-58	Kinase Assay
IKKtide	I33-58	Kinase Assay
IRS1 (Y608) Peptide	I40-58	Kinase Assay
JAK3tide	J03-58	Kinase Assay
LKBtide	L15-58	Kinase Assay
LRRKtide	L10-58	Kinase Assay
MARCKS Peptide	M02-58	Kinase Assay
METTL12 Substrate	M332-58	
Micro2 Peptide	M08-58	Kinase Assay
MLC Peptide	M09-58	Kinase Assay
Modified AKT Substrate	A05-58B	Kinase Assay
Modified AKT Substrate II	A05-58C	Kinase Assay
Modified AXLtide Substrate	A16-58B	Kinase Assay
Modified EIF2S Peptide	E23-58B	Kinase Assay
Modified PKA Substrate	C01-58B	Kinase Assay
Modified SGKtide	S08-58B	Kinase Assay

# Peptide Substrates

Product Name	Cat# No	Application
MRCL3 Peptide	M56-58	Kinase Assay
p38 Substrate	P03-58	Kinase Assay
PAKtide	P08-58	Kinase Assay
PDHKtide	P57-58	Kinase Assay
PDKtide	P10-58	Kinase Assay
PKA Substrate	C01-58	Kinase Assay
PKCtide	P15-58	Kinase Assay
PLKtide	P41-58	Kinase Assay
PLpro FRET Substrate	PL01-58	COVID-19 Assays
Poly (4:1 Glu, Tyr) Peptide	P61-58	Kinase Assay
Poly (6:2:5:1 Ala,Glu,Lys,Tyr) Peptide	P60-58	Kinase Assay
PP2B Substrate	P51-58	Phosphatase Assay
PS7-CTD Peptide	R06-58	
RS Repeat Peptide	R55-58	Kinase Assay
RSK Substrate	S06-58	Kinase Assay
S6K Substrate	S05-58	Kinase Assay
SAMStide	S07-58	Kinase Assay
SGKtide	S08-58	Kinase Assay
SRC Substrate	S30-58	Kinase Assay
TGFBR1 Peptide	T36-58	Peptide Substrate
Thr-phosphopeptide	T69-58	Phosphatase Assay
Thr-phosphopeptide-3	T72-58	Phosphatase Assay
Tyr-phosphopeptide-2	T70-58	Phosphatase Assay
Tyrosine Kinase Substrate-3	T71-58	Kinase Assay
ULKtide	U01-58	Assay
ZIPtide	Z16-58	Kinase Assay

# Protein Substrates

Product Name	Cat# No	Application
ARPP19 Protein	A76-54G	Western Blot
Casein Protein	C03-54N	Kinase Assay, Western Blot
Casein, Dephosphorylated	C03-54BN	Kinase Assay, Western Blot
DUB Substrate I	U06-57	
DUB Substrate II	S293-57	
DUB Substrate III	N263-57	
Histone H1 Protein	H10-54N	Kinase Assay, Western Blot
LC20 Protein	M89-54G	Kinase Assay, Western Blot
MBP Protein	M42-51N	Kinase Assay, Western Blot
Moesin Protein	M52-55G	Kinase Assay, Western Blot
PDHA1 Protein	P57-55H	Kinase Assay, Western Blot

**PAD Kits**



# PAD Kits

## Detection Kit

Product Name	Cat# No
P312-863	PAD Detection Kit

## Inhibitor Screening Kit

Product Name	Cat# No
P37C-863	PAD Cocktail Inhibitor Screening Kit
P318BG-863	PAD2 (human) Inhibitor Screening Kit
P310G-863	PAD1 Inhibitor Screening Kit
P310DG-863	PAD4 Inhibitor Screening Kit
P310CG-863	PAD3 Inhibitor Screening Kit
P310BG-863	PAD2 (mouse) Inhibitor Screening Kit

# CAN'T FIND THE RIGHT PRODUCT?

## SignalChem Custom Services

SignalChem understands the unique needs of scientists and researchers all over the world. To best serve those unique needs, we offer Custom Services. With our Custom Services, you can customize any product as per your requirement, expect lot-to-lot consistency and get it delivered in a competitive lead time.



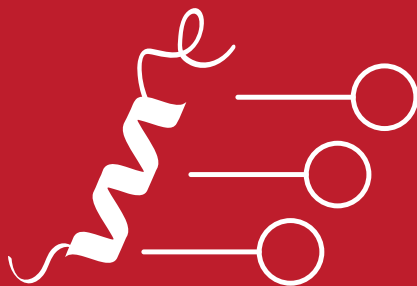
### Compound profiling

The compound selectivity profiling service offered by SignalChem is a cost-effective and convenient approach to the drug discovery continuum with only a two-week turnaround time.

### Custom Protein development

Depending upon your needs, SignalChem will work with you to custom develop the target to your desired specifications within a realistic and suitable timeframe.

Our products have been used successfully in various applications, including high-throughput screening, activity assays, binding assays, cell-based assays, protein-protein interaction studies, and structural studies.



EMAIL [info@signalchem.com](mailto:info@signalchem.com)

AND FIND THE RIGHT PRODUCT FOR YOUR RESEARCH

# SignalChem Recombinant Antibody Services

SignalChem offers recombinant antibody (rAb) productions which are generated in-vitro using cloned genes. Our custom rAbs are ethically produced without harming animals, which shows significant increase in interest and demand.

## The SignalChem Advantage

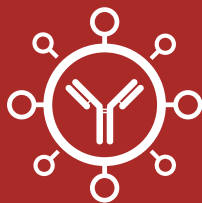
- Over 1000 in-house manufactured Active Kinases
- Comprehensive selection of Signaling Proteins, Tau Proteins, Antibodies, Compounds, and Reagents
- In-house custom services for all research products & +300 IVD products related to CRISPR, COVID -19, and Molecular Diagnostics

## Custom rAb Features

\*For all Products (unless specified)



Complete  
Expression Set



Versatility



Antibody  
Validation



Scalable  
Production

## What We Offer

- **A complete set of expression system** – Multiple expression systems - E.coli, yeast, insect, plant, and mammalian cells. Multiple Tag options - GST, HIS, MBP, TST, and others. Our expertise includes producing antibodies and their derivatives in different expression systems like mammalian cells, yeast, or E.coli.
- **Functional proteins checked for their activity, purity and antibodies** – We can also assay by Western Blot, specificity, and cross-reactivity tests, blocking function, flow cytometry staining, and other functional assays.
- **Antibody validation** – Available
- **Scalable production** – From microgram to milligram with quick turnaround time. Depending on your requirements, we can tailor the project to suit your needs.

# DISTRIBUTOR LIST

Country	Company	
All Other Countries	SignalChem Inc	
Albania	Biomedica Shpk	
Australia	Assay Matrix Pty	
Austria	Biomedica Medizinprodukte GmbH	
Belgium	DivBioScience	Bio-Connect B.V.
Bosnia and Herzegovina	Biomedica d.o.o.	
Bulgaria	Biomedica Bulgaria Ltd.	
Canada	Signalchem	Cedarlane Laboratories
China	SignalChem Biotechnologies Corporation	
Czech Republic	Biomedica CS, s.r.o	
Denmark	Nordic BioSite ApS	
Estonia	Nordic BioSite Oy	
Finland	Nordic BioSite Oy	
France	Interchim	
Germany	Biozol Diagnostica Vertrieb GmbH	
Hungary	Biomedica HungÁria Kft.	
Iceland	Nordic BioSite AB	
India	Everon Life Sciences	Biogenuix Medsystems Pvt. Ltd
Indonesia	PT Sciencewerke	
Ireland	Stratech Scientific	
Israel	Enco Scientific Services Ltd.	
Italy	DBA Italia SRL	Prodotti Gianni
Japan	Cosmo Bio	Iwai Chemicals
Latin America	ALLSCIENCE LLC	
Latvia	Nordic BioSite Oy	
Lithuania	Nordic BioSite Oy	
Luxembourg	DivBioScience	Bio-Connect B.V.
Malaysia	Sciencewerke Sdn Bhd	
Mexico	Ciencia y Biotecnologia	
Morocco	CLINI SCIENCES	
Netherlands	Sanbio B. V.	Bio-Connect B.V.
	DivBioScience	
New Zealand	Assay Matrix Pty	
North Macedonia	Biomedica dijagnostika Ltd	
Norway	Nordic BioSite AS	
Poland	Biomedica Poland Sp. z o.o.	
Portugal	Labclinics, SA	Clini Sciences
Republic of Korea	Koram Biotech	

Country	Company	
Romania	Biomedica Medizinprodukte Romania SRL	S.C. C.Y.B.E.R. Srl Romania
Serbia	Biomedica MP d.o.o.	
Singapore	Sciencewerke Pte Ltd	
Slovakia (Slovak Republic)	Biomedica Slovakia s.r.o.	
Slovenia	Biomedis M.B.	
Spain	Labclinics, SA	Condalab
Sweden	Nordic BioSite AB	
Switzerland	Lucerna Chem AG	LubioScience GmbH
Taiwan	Bio-Genesis Technologies	
Thailand	PACIFIC SCIENCE CO.,LTD.	
Tunisia	CLINI SCIENCES	
United Kingdom	Stratech Scientific	
United States	SignalChem Inc	Cedarlane Laboratories
Vietnam	Truong Bio, Inc.	

For more information, please visit [www.signalchem.com](http://www.signalchem.com)

VISIT US AT  
**SIGNALCHEM.COM**  
AND EXPERIENCE THE **POWER** OF  
**WORKING TOGETHER**



SignalChem Biotech Inc.

1.866.954.6273  
Richmond BC, Canada  
[www.signalchem.com](http://www.signalchem.com)

Last updated Aug 2023