

ACTIVE KINASE MUTANT

www.signalchem.com

Product Name	Catalog #
ABL1 Mutant (E255K), Active	A03-12BG
ABL1 Mutant (E255V), Active	A03-12HH
ABL1 Mutant (G250E), Active	A03-12CG
ABL1 Mutant (T315I), Active	A03-12DG
ABL1 Mutant (Y253F), Active	A03-12EG
ABL1 Mutant (Y253H), Active	A03-12GH
AKT1 Mutant (E17K), Active	A16-12G
AKT2 Mutant (E17K), Active	A17-12G
AKT2 Mutant (R274H), Active	A17-12BG
AKT3 Mutant (E17K), Active	A18-12G
AKT3 Mutant (G171R), Active	A18-12BG
ALK Mutant (C1156Y), Active	A19-12DG
ALK Mutant (F1174L), Active	A19-12EG
ALK Mutant (F1174S), Active	A19-12FG
ALK Mutant (G1202R), Active	A19-12HG
ALK Mutant (G1269A), Active	A19-12JG
ALK Mutant (G1269S), Active	A19-12KG
ALK Mutant (L1152R), Active	A19-12CG
ALK Mutant (L1196M), Active	A19-12GG
ALK Mutant (R1275Q), Active	A19-12LG
ALK Mutant (S1206R), Active	A19-12IG
ALK Mutant (T1151M), Active	A19-12BG
ALK Mutant (T1151_L1152insT), Active	A19-12G
ALK2 Mutant (PF197_198L), Active	A06-12FG
ALK2 Mutant (Q207D), Active	A06-13CG
ALK2 Mutant (Q207E), Active	A06-13DG
ALK2 Mutant (R206H), Active	A06-12BG
ALK3 Mutant (Q233D), Active	B04-12BG
ALK3 Mutant (R486Q), Active	B04-12CG
AXL Mutant (R499C), Active	A34-12BG
BCR-RET Mutant, Active	R02-19EG
BRAF Mutant (K601E), Active	B08-12BG
BRAF Mutant (T599_V600insT), Active	B08-13BG
BRAF Mutant (V599E), Active	B08-13G
BRAF Mutant (V600E), Active	B08-12G
BTK Mutant (C481S), Active	B10-12CG

Product Name	Catalog #
BTK Mutant (P190K), Active	B10-12BG
c-KIT Mutant (A829P), Active	C06-12HG
c-KIT Mutant (D557-558), Active	K06-12WG
c-KIT Mutant (D816E), Active	C06-12IG
c-KIT Mutant (D816F), Active	C06-12JG
c-KIT Mutant (D816H), Active	C06-12KG
c-KIT Mutant (D816V), Active	C06-12LG
c-KIT Mutant (D816Y), Active	C06-12MG
c-KIT Mutant (D820E), Active	K06-12BG
c-KIT Mutant (D820Y), Active	K06-12CG
c-KIT Mutant (K642E), Active	C06-12NG
c-KIT Mutant (L576P), Active	K06-12UG
c-KIT Mutant (N822K), Active	K06-12DG
c-KIT Mutant (T670E), Active	K06-12EG
c-KIT Mutant (T670I), Active	C06-12FG
c-KIT Mutant (V559A), Active	C06-12OG
c-KIT Mutant (V559D T670I), Active	K06-12QG
c-KIT Mutant (V559D V654A), Active	K06-12RG
c-KIT Mutant (V559D), Active	C06-12PG
c-KIT Mutant (V560G D816V), Active	C06-12TG
c-KIT Mutant (V560G N822K), Active	K06-12VG
c-KIT Mutant (V560G), Active	C06-12SG
c-KIT Mutant (V654A), Active	K06-13G
c-KIT Mutant (Y823D), Active	K06-12GG
c-MER Mutant (A708S), Active	M51-12BG
DDR1 Mutant (S496A), Active	D05-12G
DDR2 Mutant (T654M), Active	D06-12G
EGFR Mutant (d746), Active	E10-12IG
EGFR Mutant (d746-750 T790M) Protein	E10-12KG
EGFR Mutant (d746-750), Active	E10-12JG
EGFR Mutant (d747-749 A750P), Active	E10-12MG
EGFR Mutant (d747-749), Active	E10-12LG
EGFR Mutant (d747-752 P753S), Active	E10-12NG
EGFR Mutant (d752-759), Active	E10-12OG
EGFR Mutant (D761Y), Active	E10-12QG
EGFR Mutant (D770GY), Active	E10-13KG

Active Kinase Mutant

Product Name	Catalog #
EGFR Mutant (D770_N771insGF) Protein	E10-13IG
EGFR Mutant (G719C), Active	E10-12FG
EGFR Mutant (G719D), Active	E10-12GG
EGFR Mutant (G719S), Active	E10-12HG
EGFR Mutant (H773_V774insNPH) Protein	E10-13OG
EGFR Mutant (L747S), Active	E10-12PG
EGFR Mutant (L858R), Active	E10-12BG
EGFR Mutant (L861Q), Active	E10-12CG
EGFR Mutant (P772_H773insNPH), Active	E10-13MG
EGFR Mutant (T790M L858R), Active	E10-12DG
EGFR Mutant (T790M), Active	E10-12EG
EGFR Mutant (V769_D770insASV) Protein	E10-13DG
EML4-ALK Mutant, Active	A19-19G
FGFR1 Mutant (V561M), Active	F04-13G
FGFR1OP-FGFR1 Mutant, Active	F04-19BG
FGFR2 Mutant (E565G), Active	F05-12CG
FGFR2 Mutant (K526E), Active	F05-12BG
FGFR2 Mutant (K641R), Active	F05-12DG
FGFR2 Mutant (K659N), Active	F05-12EG
FGFR2 Mutant (N549H), Active	F05-12G
FGFR2 Mutant (R612T), Active	F05-12GG
FGFR2 Mutant (V564F), Active	F05-12FG
FGFR3 Mutant (G697C), Active	F06-12FG
FGFR3 Mutant (K650E), Active	F06-12CG
FGFR3 Mutant (K650M), Active	F06-12DG
FGFR3 Mutant (K650Q), Active	F06-12EG
FGFR3 Mutant (V555M), Active	F06-12GG
FGFR4 Mutant (N535K), Active	F07-12G
FGFR4 Mutant (V550E), Active	F07-12BG
FGFR4 Mutant (V550L), Active	F07-12CG
FGFR4 Mutant (V550M), Active	F07-12DG
FIP1L1-PDGFR alpha Mutant, Active	P12-19BG
FLT3 Mutant (D835Y), Active	F12-12BG
GOPC-ROS1 Mutant, Active	R14-19BG
HER2 Mutant (A775_G776insYVMA), Active	E27-13BG
HER2 Mutant (d755-759), Active	E27-12BG
HER2 Mutant (G776VC), Active	E27-13EG
HER2 Mutant (P780_Y781insGSP), Active	E27-13DG
HER2 Mutant (V777_G778insCG), Active	E27-13CG
KSR1 Mutant (A635F), Active	K07-12G
KSR1 Mutant (L639F), Active	K07-12BG
KSR2 Mutant (R676S), Active	K08-12BG
LRRK2 Mutant (D1994A), Active	L10-12FG
LRRK2 Mutant (G2019S), Active	L10-12GG
LRRK2 Mutant (G2385R), Active	L10-12IG
LRRK2 Mutant (I2020T), Active	L10-12HG
LRRK2 Mutant (R1441C), Active	L10-12G
LRRK2 Mutant (R1441G), Active	L10-12BG
LRRK2 Mutant (R1441H), Active	L10-12CG
LRRK2 Mutant (Y1699C), Active	L10-12DG
LRRK2 Mutant (Y1699G), Active	L10-12EG
MELK Mutant (T460M), Active	M50-12G

Product Name	Catalog #
MET Mutant (D1228H), Active	M52-12HG
MET Mutant (D1228N), Active	M52-12IG
MET Mutant (F1200I), Active	M52-12GG
MET Mutant (M1250T), Active	M52-12BG
MET Mutant (P991S), Active	M52-12CG
MET Mutant (T1173I), Active	M52-12FG
MET Mutant (T992I), Active	M52-12DG
MET Mutant (V1092I), Active	M52-12EG
MET Mutant (Y1230A), Active	M52-12JG
MET Mutant (Y1230C), Active	M52-12KG
MET Mutant (Y1230D), Active	M52-12LG
MET Mutant (Y1230H), Active	M52-12MG
MET Mutant (Y1235D), Active	M52-12NG
NPM1-ALK Mutant (F1174L), Active	A19-19EG
NPM1-ALK Mutant, Active	A19-19BG
PDGFR alpha Mutant (D842V), Active	P12-12BG
PDGFR alpha Mutant (T674I), Active	P12-12CG
PI3K Mutant (p110 alpha(E542K)/p85 alpha), Active	P27-12DH
PI3K Mutant (p110 alpha(E545K)/p85 alpha), Active	P27-15H
PI3K Mutant (p110alpha(N345K)/p85alpha), Active	P27-12CH
PKD2 Mutant (G848E), Active	P76-12BG
PKD2 Mutant (G870E), Active	P76-12CG
PTC1 Mutant (CCDC6-RET), Active	R02-19BG
PTC2 Mutant (PRKAR1A-RET), Active	R02-19CG
PTC3 Mutant (NCOA4-RET), Active	R02-19DG
RET Mutant (A883F), Active	R02-12MG
RET Mutant (E762Q), Active	R02-12FG
RET Mutant (G691S), Active	R02-12DG
RET Mutant (L790F), Active	R02-12LG
RET Mutant (M918T), Active	R02-12JG
RET Mutant (R749T), Active	R02-12EG
RET Mutant (R813Q), Active	R02-12HG
RET Mutant (S891A), Active	R02-12IG
RET Mutant (V778I), Active	R02-12KG
RET Mutant (V804L), Active	R02-12BG
RET Mutant (V804M), Active	R02-12GG
RET Mutant (Y791F), Active	R02-12CG
RSK2 Mutant (I416V), Active	R17-12BG
RSK2 Mutant (L608F), Active	R17-12CG
RSK2 Mutant (Y483C), Active	R17-12DG
TGFBR1 Mutant (T204D), Active	T07-13G
TIE2 Mutant (A1124V), Active	T04-12GG
TIE2 Mutant (P883A), Active	T04-12CG
TIE2 Mutant (R849W), Active	T04-12BG
TIE2 Mutant (Y1108F), Active	T04-12FG
TIE2 Mutant (Y897C), Active	T04-12DG
TIE2 Mutant (Y897S), Active	T04-12EG
TPM3-ALK Mutant, Active	A19-19DG
TPM3-ROS1 Mutant, Active	R14-19CG
TPM3-TRKA Mutant, Active	N16-19DG
TPR-TRKA Mutant (TRK-T1), Active	N16-19BG
ZAP70 Mutant (Y319F), Active	Z02-12BG