



# NEUROBIOLOGY

Product Name	Catalog #
14-3-3 alpha/beta Protein	Y71-30G
14-3-3 theta Protein	Y84-30G
ABL1, Active	A03-11H
ABL2, Active	A04-11H
ACK, Active	A05-11G
AKT1, Active	A16-10G
AKT2, Active	A17-10H
AKT3, Active	A18-10G
ALK, Active	A19-11G
ALK1, Active	A09-11G
ALK2, Active	A06-11G
ALK3 (BMPR1A), Active	B04-11G
ALK4, Active	A07-11G
ALK6 (BMPR1B), Active	B05-11G
ATG13 Protein	A113-30G
BID Protein	B59-30G
c-JUN Protein	J05-55G
c-KIT, Active	K06-11G
CAMK1 beta, Active	C08-10G
CAMK1 delta, Active	C09-10G
CAMK2 beta, Active	C12-10H
CAMK2 gamma, Active	C14-10G
CAMKK1, Active	C17-10G
CAMKK2, Active	C18-10G
CaMKPase Protein	P06-34G
CASK, Active	C19-11G
CK1 alpha 1, Active	C64-10G
Cofilin 1 Protein	C37-54H
CREB1 Protein	C49-55G
CRYAA Protein	C256-30H
DAPK1, Active	D01-11G
DAPK2, Active	D02-10G
DCAMKL1, Active	D14-10G
DCAMKL2, Active	D15-11G
DNMT1 Protein	D351-31G
DNMT3L, Active	D353-380CG
DUSP22 (MKPX), Active	D22-20G
DYRK2, Active	D10-10G
DYRK3, Active	D11-11G
EEF2K, Active	E01-10G
EHMT2, Active	E318-381G
EIF2AK1 (HRI), Active	H07-10G
EIF2AK4 (GCN2), Active	E12-11G
EPHA1, Active	E13-11G
EPHA2, Active	E14-11G
EPHA4, Active	E16-11G
EPHB1, Active	E21-11G
EPHB2, Active	E22-11G
EPHB4, Active	E24-11G

Product Name	Catalog #
ERK1, Active	M29-10G
ERK2, Active	M28-10G
FER, Active	F16-11G
FGFR1 (FLT2), Active	F04-11G
FYN A, Active	F15-10G
GAB1 Protein	G03-30H
GRB7 Protein	G45-30GG
GRK1, Active	R03-10G
GRK3, Active	A15-11G
GSK3 alpha, Active	G08-10G
GSK3 beta, Active	G09-10G
HGK, Active	M26-11G
HO1 Protein	H20-30G
IKK alpha, Active	C51-10G
IKK beta, Active	I03-10BG
JNK1, Active	M33-10G
JNK2, Active	M34-10BG
KAT2B (PCAF), Active	K311-381BG
KSR2, Active	K08-11G
LRRK2, Active	L10-11G
MAPKAPK2, Active	M40-11G
MAPKAPK2, Active	M40-11H
MARK1, Active	M43-10G
MARK2, Active	M44-10G
MARK3, Active	M45-10G
MARK4, Active	M46-10G
MBP Protein	M42-54G
MBP Protein	M42-51N
MEK1, Active	M02-10G
MEK1 (EE), Active	M02-12G
MEK2, Active	M03-10G
MELK, Active	M50-11G
MINK1, Active	M53-11G
MSK1, Active	R19-10G
MST1, Active	S25-10G
MUSK, Active	M60-11G
MYO3 beta, Active	M66-11G
N6AMT2 Protein	N355-30BH
NIK, Active	M22-11G
p38 alpha, Active	M39-10BG
p53 Protein	P05-30BG
p63 Protein	P06-30G
PAK1, Active	P02-10BG
PASK, Active	P08-11G
PDK1, Active	P14-10H
PI3K (p110 alpha/p85 alpha), Active	P27-10H
PI3K (p110 alpha/p85 alpha), Active	P27-18H
PI3K (p110 delta/p85 alpha), Active	P30-10H
PI3K (p120 gamma), Active	P29-10H

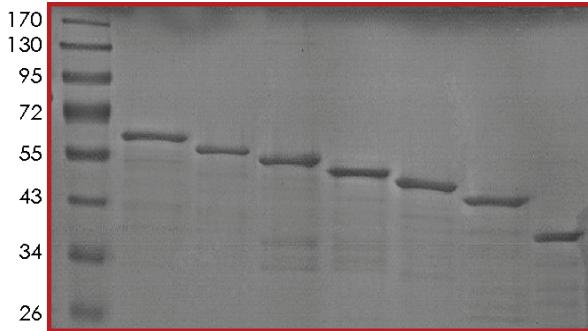
Product Name	Catalog #
PI4K2A, Active	P21-10G
PI4K2B, Active	P22-10G
PI4KB, Active	P32-10G
PIK3C2A, Active	P23-11G
PIK3C2G, Active	P25-11G
PIK3C3 (Vps34), Active	P26-10G
PIM1, Active	P35-10G
PIM3, Active	P37-10G
PIP4K2A, Active	P76-10AG
PIP4K2C, Active	P76-10CG
PIP5K1B, Active	P16-10BG
PIP5K1C, Active	P16-10CG
PIP5K3 Protein	P17-31G
PKAc alpha, Active	P51-10G
PKD2, Active	P76-10G
PKN2/PRK2, Active	P71-10G
PKN3/PRK3, Active	P69-10G
PRKG1, Active	P78-10BG
PRMT1, Active	P365-380G
PRMT3, Active	P365-380CG
PSKH1 Protein	P89-34G
PTEN, Active	P23-20G
PTPN7 (LC-PTP), Active	P34-20G
PTPRA (LRP), Active	P48-21G
PYK2, Active	P92-11H
RAC2 Protein	R04-30H
RAF1 (EE), Active	R01-13G
RAS Protein	R05-30G
RET, Active	R02-11G
RhoD Protein	R44-30DH
ROR1, Active	R12-11G
ROR2, Active	R13-11G
RSK1, Active	R15-10G
SETD7, Active	S344-380G
SETD8 Protein	S345-30G
SGK1, Active	S06-11G
SOK1, Active	S43-10G
SPHK1, Active	S17-10G
STK9 Protein	C46-31G
STK32C (YANK3), Active	Y05-11G
STK33, Active	S46-10G
STK36, Active	S48-11G
STK39 (STLK3), Active	S51-10H
SUV39H1 Protein	S350-30G
TIE 2, Active	T04-11G
TLK1, Active	T10-10G
TNIK, Active	T27-11G
TYRO3, Active	T22-11G
XIAP Protein	X08-30G

# TAU Proteins

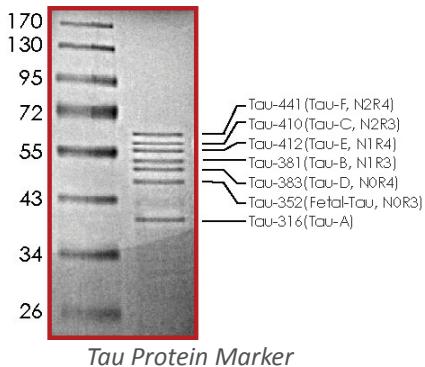
It is well established that many neurodegenerative diseases result from the selective loss of neurons within central nervous system; however, the precise mechanism(s) leading to this susceptibility remains somewhat elusive. In a portion of neurological disorders, biochemical alterations in a single protein are associated with the observed clinical manifestations.

Tau is one such protein that is linked to the pathogenesis of a number of neurodegenerative diseases. Tau is a microtubule-associated protein that is heat resistant. Under normal circumstances, this phosphoprotein plays a role in regulating microtubule polymerization and stabilization. In neurodegenerative diseases, the Tau protein is known to be either mutated or possess amino acid substitutions; to date over 30 distinct mutations have been noted. These genetic modifications in Tau likely alter its conformation and solubility leading to its aggregation and the formation of fibrils.

## Highly Quality Tau Proteins Available for Research



The image of seven Tau proteins, from left to right: Tau-441, Tau-410, Tau-412, Tau-381, Tau-383, Tau-352 and Tau-316.



Tau Protein Marker

Product Name	Catalog #	Product Name	Catalog #	Product Name	Catalog #
Tau Protein Marker	T08-07N	Tau-441 (S198E) Protein	T08-53N	Tau-441 (L266V) Protein	T08-56CN
Tau-316 Protein	T02-54N	Tau-441 (1-421) Protein	T08-55BN	Tau-441 (G272V) Protein	T08-56DN
Tau-352 Protein	T03-54N	Tau-441 (1-391) Protein	T08-55CN	Tau-441 (N279K) Protein	T08-56EN
Tau-381 Protein	T04-54N	Tau-441 (151-441) Protein	T08-55DN	Tau-441 (P301L) Protein	T08-56FN
Tau-383 Protein	T05-54N	Tau-441 (151-421) Protein	T08-55EN	Tau-441 (P301S) Protein	T08-56GN
Tau-383 (N352H) Protein	T05-56N	Tau-441 (151-391) Protein	T08-55FN	Tau-441 (S305N) Protein	T08-56HN
Tau-410 Protein	T06-54N	Tau-441 (231-441) Protein	T08-55GN	Tau-441 (S352L) Protein	T08-56KN
Tau-412 Protein	T07-54N	Tau-441 (231-421) Protein	T08-55HN	Tau-441 (R406W) Protein	T08-56MN
Tau-441 Protein	T08-54H	Tau-441 (231-391) Protein	T08-55IN	Tau-441 (K257T) Protein	T08-56N
Tau-441 Protein	T08-54N	Tau-441 (99-441) Protein	T08-55JN	Tau-441, BRSK1-phosphorylated	T08-50N
Tau-441 (dN296) Protein	T08-52BN	Tau-441 (127-421) Protein	T08-55KN	Tau-441, BRSK2-phosphorylated	T08-50BN
Tau-441 (dK280) Protein	T08-52N	Tau-441 (216-391) Protein	T08-55LN	Tau-441, CAMK2-phosphorylated	T08-50CN
Tau-441 (S198A) Protein	T08-53BN	Tau-441 (50-441) Protein	T08-55MN	Tau-441, GSK3beta-phosphorylated	T08-50FN
Tau-441 (S199E) Protein	T08-53CN	Tau-441 (244-372) Protein	T08-55N	Tau-441, PHKG2-phosphorylated	T08-50KN
Tau-441 (S214A) Protein	T08-53DN	Tau-441 (50-421) Protein	T08-55NN	Tau-441, PKA-phosphorylated	T08-50LN
Tau-441 (S404A) Protein	T08-53HN	Tau-441 (50-391) Protein	T08-55ON	Tau-441, TTBK1-phosphorylated	T08-50ON