



# SER/THR KINASES

www.signalchem.com

Product Name	Cat #	Product Name	Cat #	Product Name	Cat #
ACVR2A Protein	A35-35H	CAMKK1, Active	C17-18G	DRAK1 (STK17A), Active	S33-10G
AKT1, Active	A16-10G	CAMKK2, Active	C18-10G	DRAK2 (STK17B), Active	S34-10G
AKT1 Mutant (E17K), Active	A16-12G	CASK, Active	C19-11G	DYRK2, Active	D10-10G
AKT2, Active	A17-10G	CDC7 Protein	C26-30G	DYRK3, Active	D11-11G
AKT2, Active	A17-10H	CDC7/DBF4, Active	C26-10G	EEF2K, Active	E01-10G
AKT2 Mutant (E17K), Active	A17-12G	CDC25A, Active	C04-20G	EIF2AK1 (HRI), Active	H07-10G
AKT2 Mutant (R274H), Active	A17-12BG	CDC25B, Active	C04-20BG	EIF2AK2, Active	P80-11G
AKT3, Active	A18-10G	CDC25C, Active	C04-20CG	EIF2AK3, Active	E11-11G
AKT3 Mutant (G171R), Active	A18-12BG	CDK1/CyclinA1, Active	C22-18BG	EIF2AK4 (GCN2), Active	E12-11G
ALK, Active	A19-11G	CDK1/CyclinA2, Active	C22-18G	ERK1, Active	M29-10G
ALK Mutant (F1174S), Active	A19-12FG	CDK1/CyclinB1, Active	C22-10G	ERK1, Active	M29-10U
ALK Mutant (L1196M), Active	A19-12GG	CDK2/CyclinA1, Active	C29-10BG	ERK2, Active	M28-10G
ALK Mutant (S1206R), Active	A19-12IG	CDK2/CyclinA2, Active	C29-10G	ERK5, Active	M32-10G
ALK1, Active	A09-11G	CDK2/CyclinE1, Active	C29-18G	FASTK, Active	F01-10G
ALK2, Active	A06-11G	CDK2/CyclinO, Active	C29-19G	GCK, Active	M24-10G
ALK2 Mutant (R206H), Active	A06-12BG	CDK2 Protein	C29-30G	GLK, Active	M25-11G
ALK3 (BMPR1A), Active	B04-11G	CDK3/CyclinE1, Active	C30-10G	GRK1, Active	R03-10G
ALK4, Active	A07-11G	CDK4/CyclinD1, Active	C31-10G	GRK2, Active	A14-10G
ALK6 (BMPR1B), Active	B05-11G	CDK4/CyclinD3, Active	C31-18G	GRK3, Active	A15-11G
AMPK (A1/B1/G1), Active	P47-10H	CDK5/p25, Active	C33-10G	GRK5, Active	G03-10G
AMPK (A1/B1/G2), Active	P55-10H	CDK5/p35, Active	C33-10BG	GRK6, Active	G04-10G
AMPK (A1/B1/G3), Active	P56-10H	CDK6/CyclinD1, Active	C35-18H	GRK7, Active	G05-10G
AMPK (A1/B2/G1), Active	P50-10H	CDK6/CyclinD3, Active	C35-10H	GSK3 alpha, Active	G08-10G
AMPK (A1/B2/G2), Active	P45-10G	CDK7/CyclinH1/MNAT1, Active	C36-10H	GSK3 beta, Active	G09-10G
AMPK (A1/B2/G3), Active	P83-10G	CDK9/CyclinK, Active	C40-10G	GSK3 beta, Active	G09-10H
AMPK (A2/B1/G1), Active	P48-10H	CHK1, Active	C47-10G	Haspin (GSG2), Active	G10-11G
AMPK (A2/B1/G2), Active	P84-10GH	CHK1, Active	C47-10H	HGK, Active	M26-11G
AMPK (A2/B1/G3), Active	P85-10GH	CHK2, Active	C48-10G	HIPK1, Active	H03-11G
AMPK (A2/B2/G1), Active	P49-10H	CK1 alpha 1, Active	C64-10G	HIPK2, Active	H04-11G
AMPK (A2/B2/G2), Active	P54-10H	CK1 delta, Active	C65-10G	HIPK3, Active	H05-13H
AMPK (A2/B2/G3), Active	P46-10GH	CK1 epsilon, Active	C66-10G	HIPK4, Active	H06-10G
BRAF, Active	B08-11G	CK1 gamma 1, Active	C68-11G	ICK, Active	I01-10G
BRSK1, Active	B13-10G	CK1 gamma 2, Active	C68-10BG	IKK alpha, Active	C51-10G
BRSK2, Active	B14-10G	CK1 gamma 3, Active	C68-10CG	IKK beta, Active	I03-10BG
BUB1B, Active	B12-11G	CK2 alpha 1, Active	C70-10G	IRAK2, Active	I10-10BG
CAMK1, Active	C07-10G	CK2 alpha 2, Active	C71-10G	IRAK4, Active	I12-10G
CAMK1 beta, Active	C08-10G	CLK1, Active	C57-11G	JNK1, Active	M33-10G
CAMK1 delta, Active	C09-10G	CLK2, Active	C58-11G	JNK2, Active	M34-10BG
CAMK1 gamma, Active	C10-10BG	CLK3, Active	C60-10G	JNK3, Active	M35-10BG
CAMK1 gamma, Active	C10-11G	COT, Active	M16-11G	KHS1, Active	M27-10G
CAMK2 alpha, Active	C11-10G	DAPK1, Active	D01-11G	KSR1, Active	K07-11G
CAMK2 beta, Active	C12-10H	DAPK2, Active	D02-10G	KSR2, Active	K08-11G
CAMK2 delta, Active	C13-10G	DAPK3, Active	D03-10G	KSR2 Mutant (R676S), Active	K08-12BG
CAMK2 gamma, Active	C14-10G	DCAMK1, Active	D14-10G	LIMK1, Active	L04-11G
CAMK4, Active	C15-10G	DCAMK2, Active	D15-11G	LOK, Active	S29-11G
CAMKK1, Active	C17-10G	DMPK, Active	D07-10G	LTK, Active	L11-11G

# SER/THR Kinases

Product Name	Cat #
MAPKAPK2, Active	M40-11G
MAPKAPK2, Active	M40-11H
MAPKAPK3, Active	M41-10G
MAPKAPK5, Active	M42-10G
MARK1, Active	M43-10G
MARK2, Active	M44-10G
MARK3, Active	M45-10G
MARK4, Active	M46-10G
MEK1, Active	M02-10G
MEK1 (EE), Active	M02-12G
MEK2, Active	M03-10G
MEKK1, Active	M09-11G
MEKK2, Active	M10-10G
MEKK3, Active	M11-10G
MEKK6, Active	M14-11G
MELK, Active	M50-11G
MELK Mutant (T460M), Active	M50-12G
MINK1, Active	M53-11G
MLCK, Active	M62-11G
MLK1, Active	M17-11G
MLK2, Active	M18-11G
MLK3, Active	M19-11G
MLK4, Active	M48-11G
MNK1, Active	M54-10G
MNK2, Active	M55-10G
MRCK alpha, Active	C27-11G
MRCK beta, Active	C28-11G
MSK1, Active	R19-10G
MSK2, Active	R18-10G
MSSK1, Active	S41-10G
MYLK2, Active	M63-10H
MYLK3, Active	M64-10G
MYO3 beta, Active	M66-11G
NDR, Active	S49-10G
NDR2, Active	S50-10G
NEK1, Active	N02-11G
NEK2, Active	N03-10G
NEK3, Active	N05-10G
NEK5, Active	N07-11G
NEK6, Active	N08-10G
NEK7, Active	N09-10G
NIK, Active	M22-11G
NIM1, Active	N14-10G
NLK, Active	N15-10G
NUAK1, Active	N19-10G
NUAK2, Active	N20-10G
p38 alpha, Active	M39-10BG
p38 beta, Active	M36-10BG
p38 delta, Active	M38-10BG
p38 gamma, Active	M37-10BG
p70S6K, Active	R21-10H
p70S6Kb, Active	R22-10G
PAK1, Active	P02-10BG
PAK1/CDC42, Active	P02-10G
PAK3, Active	P04-10G

Product Name	Cat #
PAK4, Active	P05-10BG
PAK6, Active	P06-10G
PAK7, Active	P07-10G
PASK, Active	P08-11G
PCTK1 (CDK16)/CyclinY, Active	P09-11G
PCTK3/CyclinY, Active	P11-10G
PDHK1, Active	P57-10G
PDHK2, Active	P58-10G
PDHK3, Active	P59-10G
PDHK4, Active	P60-10G
PDK1, Active	P14-10H
PFTK1 (CDK14)/CyclinY, Active	P15-10G
PHKG1, Active	P19-10G
PHKG2, Active	P20-10G
PIM1, Active	P35-10G
PIM2, Active	P36-10G
PIM3, Active	P37-10G
PKAc alpha, Active	P51-10G
PKAc beta, Active	P52-10G
PKAc gamma, Active	P53-10G
PKC alpha, Active	P61-10G
PKC alpha, Active	P61-18G
PKC beta I, Active	P62-10G
PKC beta II, Active	P63-10G
PKC delta, Active	P64-10G
PKC epsilon, Active	P65-10G
PKC eta, Active	P67-10G
PKC gamma, Active	P66-10G
PKC iota, Active	P68-10G
PKC mu, Active	P72-10G
PKC nu, Active	P73-10G
PKC theta, Active	P74-10G
PKC zeta, Active	P75-10G
PKD2, Active	P76-10G
PKD2 Mutant (G848E), Active	P76-12BG
PKD2 Mutant (G870E), Active	P76-12CG
PKN2/PRK2, Active	P71-10G
PKN3/PRK3, Active	P69-10G
PLK1, Active	P41-10H
PLK2, Active	P42-10G
PLK3, Active	P43-13G
PLK4, Active	P44-11G
PP1A, Active	P13-20G
PP1B, Active	P14-20G
PP2Aalpha, Active	P16-20G
PP2Calpha, Active	P02-20G
PP2Cgamma, Active	P07-20G
PRKG1, Active	P78-10BG
PRKG1, Active	P78-11G
PRKG2, Active	P79-10G
PRKX, Active	P81-10G
QIK, Active	S15-10G
RAF1 (EE), Active	R01-13G
RIPK2, Active	R08-11G
RIPK3, Active	R09-10G

Product Name	Cat #
RIPK5, Active	R27-10G
ROCK1, Active	R10-11G
ROCK2, Active	R11-11G
ROCK2, Active	R11-11H
RSK1, Active	R15-10G
RSK2, Active	R17-10G
RSK2 Mutant (I416V), Active	R17-12BG
RSK2 Mutant (Y483C), Active	R17-12DG
RSK2 Mutant (L608F), Active	R17-12CG
RSK3, Active	R16-10G
RSK4, Active	R20-10G
SGK1, Active	S06-11G
SGK2, Active	S07-10G
SGK3, Active	S08-10G
SIK, Active	S14-11H
SIK3, Active	S12-11G
SLK, Active	S11-10G
SOK1, Active	S43-10G
SRPK1, Active	S21-10G
SRPK2, Active	S22-10G
STK3, Active	S24-10G
STK19, Active	S35-10G
STK21 (CIT), Active	C52-11G
STK32B (YANK2), Active	Y04-10G
STK32C (YANK3), Active	Y05-11G
STK33, Active	S46-10G
STK36, Active	S48-11G
STK39 (STLK3), Active	S51-10H
TAK1-TAB1, Active	M15-13G
TAOK1, Active	T24-11G
TAOK2, Active	T25-11G
TAOK3, Active	T26-11G
TBK1, Active	T02-10G
TESK1 Protein	T05-34G
TESK2, Active	T06-10G
TGFBR1 (ALK5), Active	T07-11G
TGFBR2, Active	T08-11G
TLK2, Active	T11-11G
TNIK, Active	T27-11G
TOPK, Active	T14-10G
TSSK1B, Active	S40-10BG
TSSK2, Active	S38-10G
TSSK3 Protein	S39-34G
TTBK1, Active	T17-11G
TTBK2, Active	T18-11G
TTK, Active	T20-10G
ULK1, Active	U01-11G
ULK2, Active	U02-11G
ULK3, Active	U03-10H
VRK1, Active	V01-10G
VRK2, Active	V02-11G
WNK1, Active	W02-11G
ZAK, Active	Z01-10G