

TYROSINE KINASES

Receptor Tyrosine Kinases

Product Name	Cat #
ACK, Active	A05-11G
AXL, Active	A34-11H
AXL Mutant (R499C), Active	A34-12BG
c-MER, Active	M51-11G
CRKL Protein	C53-30H
DDR1, Active	D05-11G
DDR2, Active	D06-11BG
DDR2 Mutant (T654M), Active	D06-12G
EGFR, Active	E10-11G
EGFR Mutant (T790M), Active	E10-12EG
EGFR Mutant (T790M L858R), Active	E10-12DG
EGFR Mutant (L858R), Active	E10-12BG
EGFR Mutant (L861Q), Active	E10-12CG
EPHA1, Active	E13-11G
EPHA1, Active	E13-18G
EPHA2, Active	E14-11G
EPHA3, Active	E15-11G
EPHA4, Active	E16-11G
EPHA6, Active	E18-11G
EPHA7, Active	E19-11G
EPHA8 Protein	E20-35G
EPHB1, Active	E21-11G
EPHB2, Active	E22-11G
EPHB3, Active	E23-11G
EPHB4, Active	E24-11G
EPHB4, Active	E24-11H
FGFR1 (FLT2), Active	F04-11G
FGFR1 Mutant (V561M), Active	F04-13G
FGFR2, Active	F05-11G
FGFR2 Mutant (K526E), Active	F05-12BG
FGFR2 Mutant (N549H), Active	F05-12G

Product Name	Cat #
FGFR2 Mutant (E565G), Active	F05-12CG
FGFR2 Mutant (K641R), Active	F05-12DG
FGFR2 Mutant (K659N), Active	F05-12EG
FGFR3, Active	F06-11G
FGFR3 Mutant (K650E), Active	F06-12CG
FGFR3 Mutant (K650M), Active	F06-12DG
FGFR3 Mutant (K650Q), Active	F06-12EG
FGFR3 Mutant (G697C), Active	F06-12FG
FGFR4, Active	F07-11G
FGFR4 Mutant (N535K), Active	F07-12G
FGFR4 Mutant (V550E), Active	F07-12BG
FGFR4 Mutant (V550L), Active	F07-12CG
FGFR4 Mutant (V550M), Active	F07-12DG
FLT1, Active	F11-11G
FLT3, Active	F12-11G
FLT3 Mutant (D835Y), Active	F12-12BG
FLT4, Active	F13-11G
HER2, Active	E27-11G
HER4, Active	E29-11G
IGF1R, Active	I02-11H
InsR, Active	I08-11G
IRR, Active	I07-11G
KDR, Active	K01-11G
MET, Active	M52-11G
MET, Active	M52-18G
MET Mutant (P991S), Active	M52-12CG
MET Mutant (T992I), Active	M52-12DG
MET Mutant (V1092I), Active	M52-12EG
MET Mutant (T1173I), Active	M52-12FG
MET Mutant (F1200I), Active	M52-12GG
MET Mutant (D1228H), Active	M52-12HG

Product Name	Cat #
MET Mutant (D1228N), Active	M52-12IG
MET Mutant (Y1230C), Active	M52-12KG
MET Mutant (Y1230D), Active	M52-12LG
MET Mutant (Y1230H), Active	M52-12MG
MET Mutant (Y1235D), Active	M52-12NG
MET Mutant (M1250T), Active	M52-12BG
MST1, Active	S25-10G
MST3, Active	S42-11G
MST4, Active	M59-10G
MUSK, Active	M60-11G
PDGFR alpha, Active	P12-11G
PDGFR alpha, Active	P12-18G
PDGFR alpha Mutant (T674I), Active	P12-12CG
PDGFR alpha Mutant (D842V), Active	P12-12BG
PDGFR beta, Active	P13-11G
RET, Active	R02-11G
RET Mutant (Y791F), Active	R02-12CG
RET Mutant (V804L), Active	R02-12BG
RON, Active	M58-11G
ROR1, Active	R12-11G
ROR2, Active	R13-11G
ROS1, Active	R14-11G
RYK Protein	R28-31G
TEC, Active	T03-10G
TIE 2, Active	T04-11G
TLK1, Active	T10-10G
TRKA, Active	N16-11G
TRKB, Active	N17-11G
TRKC, Active	N18-11H
TYRO3, Active	T22-11G

Cytoplasmic Tyrosine Kinases

Product Name	Cat #
AATK Protein	A02-35G
ABL1, Active	A03-11H
ABL1, Active	A03-18H
ABL1 Mutant (E255K), Active	A03-12BG
ABL1 Mutant (G250E), Active	A03-12CG
ABL1 Mutant (T315I), Active	A03-12DG
ABL1 Mutant (Y253F), Active	A03-12EG
ABL2, Active	A04-11H
BLK, Active	B02-10G
BMX, Active	B07-10G
BRK, Active	P94-10G
BTK, Active	B10-10H
c-KIT, Active	K06-11BG
c-KIT, Active	K06-11G
c-KIT Mutant (V559A), Active	C06-12OG
c-KIT Mutant (V559D), Active	C06-12PG
c-KIT Mutant (V559D T670I), Active	C06-12QG
c-KIT Mutant (V559D T670I), Active	K06-12QG
c-KIT Mutant (V559D V654A), Active	C06-12RG
c-KIT Mutant (V559D V654A), Active	K06-12RG
c-KIT Mutant (V560G D816V), Active	C06-12TG
c-KIT Mutant (V560G), Active	C06-12SG
c-KIT Mutant (K642E), Active	C06-12NG
c-KIT Mutant (V654A), Active	K06-13G
c-KIT Mutant (T670E), Active	C06-12EG
c-KIT Mutant (T670E), Active	K06-12EG

Product Name	Cat #
c-KIT Mutant (T670I), Active	C06-12FG
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	C06-12BG
c-KIT Mutant (D820E), Active	K06-12BG
c-KIT Mutant (N822K), Active	C06-12DG
c-KIT Mutant (N822K), Active	K06-12DG
c-KIT Mutant (Y823D), Active	C06-12GG
c-KIT Mutant (Y823D), Active	K06-12GG
c-KIT Mutant (A829P), Active	C06-12HG
CSK, Active	C63-10G
CSK, Active	C63-10U
FAK, Active	P91-10G
FAK, Active	P91-11H
FER, Active	F16-11G
FES, Active	F03-10G
FGR, Active	F10-10G
FMS, Active	C74-11G
FRK, Active	F14-11G
FYN A, Active	F15-10G
FYN C, Unactive	F15-14G
HCK, Active	H02-11G
HERC4 Protein	H265-31G

Product Name	Cat #
HPK1, Active	M23-11G
ITK, Active	I13-11G
JAK2, Active*	J02-11G
JAK2, Active*	J02-11H
JAK3, Active	J03-11G
LCK, Active	L03-10G
LYN A, Active	L13-18G
LYN B, Active	L13-10G
PEAK1, Active	P18-11G
PEAK1, Unactive	P18-14G
PYK2, Active	P92-11H
SRC, Active	S19-10G
SRC, Active	S19-18G
SRMS, Active	S20-11G
STUB1 (CHIP) Protein	C268-30H
STYK1 (NOK) Protein	S53-35G
SYK, Active	S52-10G
TTK, Active	T20-10G
TXK, Active	T19-11G
TYK2 (JTK1), Active	T21-11G
UBA1 (UBE1) Protein	U201-30G
VLK Protein	V05-31G
WEE1 Protein	W01-35G
YES1, Active	Y01-10G
ZAP70, Active	Z02-10G

Products for All Seven Kinase Families

TK, TKL, STE, CK1, AGC, CAMK, and CMGC

⊕ **HIGH SPECIFIC ACTIVITY**
Good Signal to Noise Ratio = Greater Utility

⊕ **HIGH PURITY**
Low Non-Specific Interference = Larger Success

* This product shall not be used to commercially screen drug molecules being developed as JAK1 or JAK2 inhibitors nor for any in vitro diagnostic purposes. Any such activity will be outside the scope of the research use only label license. The product is protected by the following patents: US5,910,426; US5,852,184; US5,821,069; US5,716,818; US5,658,791.

20140115

email: ORDERS@SIGNALCHEM.COM

toll free: 1.866.9.KINASE

phone: 1.604.232.4600