

KinaseSeeker™ Full-Length Panel (112 Wild-Type Kinases)

Kinase	Group
ABL1 full-length	TK
ACVR1 intracellular module	TKL
AKT1 full-length	AGC
AKT2 full-length	AGC
AKT3 full-length	AGC
AMPKa1 full-length	CAMK
BLK full-length	TK
BTK full-length	TK
CAMK1D full-length	CAMK
CAMK1G full-length	CAMK
CAMK2A full-length	CAMK
CAMK2B full-length	CAMK
CAMK2D full-length	CAMK
CAMK2G full-length	CAMK
CAMKK1 full-length	Other
CAMKK2 full-length	Other
CASK full-length	CAMK
CDKL5 full-length	CMGC
CK1d full-length	CK1
CLK1 full-length	CMGC
CLK2 full-length	CMGC
CLK3 full-length	CMGC
CSF1R intracellular module	TK
CSK full-length	TK
DAPK1 full-length	CAMK

Kinase	Group
DDR1 intracellular module	TK
DDR2 intracellular module	TK
EGFR intracellular module	TK
EPHA1 intracellular module	TK
EPHA2 intracellular module	TK
EPHA3 intracellular module	TK
EPHA4 intracellular module	TK
EPHA5 intracellular module	TK
EPHA6 intracellular module	TK
EPHA7 intracellular module	TK
EPHA8 intracellular module	TK
EPHB3 intracellular module	TK
EPHB4 intracellular module	TK
ERBB2 intracellular module	TK
ERBB4 intracellular module	TK
FAK full-length	TK
FGFR2 intracellular module	TK
FGFR3 intracellular module	TK
FGR full-length	TK
FLT1 intracellular module	TK
FLT2 intracellular module	TK
FLT4 intracellular module	TK
FRK full-length	TK
FYN full-length	TK
GRK7 full-length	AGC

Legend: Full-Length: Construct contains Full-length kinase

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

KinaseSeeker™ Full-Length Panel (112 Wild-Type Kinases)

Kinase	Group
GSK3A full-length	CMGC
HCK full-length isoform1	TK
HCK full-length isoform2	TK
HIPK4 full-length	CMGC
IGF1R intracellular module	TK
IKKe full-length	Other
INSR intracellular module	TK
IRAK3 full-length	TKL
KIT intracellular module	TK
LCK full-length	TK
LYN full-length	TK
MAP4K2 full-length	STE
MAP4K5 full-length	STE
MAPK9 full-length	CMGC
MARK1 full-length	CAMK
MARK2 full-length	CAMK
MET intracellular module	TK
MSK1 full-length	AGC
MSK2 full-length	AGC
MST1 full-length	STE
MST2/STK3 full-length	STE
MST3 full-length	STE
MUSK intracellular module	TK
MYLK2 full-length	CAMK
MYLK4 full-length	CAMK

Kinase	Group
MYT1 full-length	Other
NTRK2 intracellular module	TK
NTRK3 intracellular module	TK
NUAK1 full-length	CAMK
PDGFRA intracellular module	TK
PDGFRB intracellular module	TK
PDK1 full-length	AGC
PFTAIRE1 full-length	CMGC
PKG1 full-length	AGC
PLK4 full-length	Other
PTK6 full-length	TK
PYK2 full-length	TK
RET intracellular module	TK
RIPK2 full-length	TKL
ROS1 intracellular module	TK
RSK1 full-length	AGC
RSK2 full-length	AGC
RSK3 full-length	AGC
RSK4 full-length	AGC
SIK1 full-length	CAMK
SIK2 full-length	CAMK
SLK full-length	STE
SNARK full-length	CAMK
SRC full-length	TK
SRC (Y530F) full-length	TK

Legend: Full-Length: Construct contains Full-length kinase

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

KinaseSeeker™ Full-Length Panel (112 Wild-Type Kinases)

Kinase	Group
STK10 full-length	STE
STK33 full-length	CAMK
TAOK2 full-length	STE
TBK1 full-length	Other
TESK2 full-length	TKL
TIE1 intracellular module	TK
TIE2 intracellular module	TK
TNK1 full-length	TK
TYRO3 intracellular module	TK
ULK2 full-length	Other
VEGFR2 intracellular module	TK
YES1 full-length	TK

To place an order or request a quote,
please contact us at:

Phone: (520) 495-0161
Email: info@luceome.com

Legend: Full-Length: Construct contains Full-length kinase
Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases