

KinaseSeeker™ TK Panel (126 Wild-Type and Full-Length Kinases)

Kinase	Group
ABL1	TK
ABL1 full-length	TK
ABL2	TK
ALK	TK
AXL	TK
BLK	TK
BLK full-length	TK
BMX	TK
BTK	TK
BTK full-length	TK
CSF1R	TK
CSF1R intracellular module	TK
CSK	TK
CSK full-length	TK
DDR1	TK
DDR1 intracellular module	TK
DDR2	TK
DDR2 intracellular module	TK
EGFR	TK
EGFR intracellular module	TK
EPHA1	TK
EPHA1 intracellular module	TK
EPHA2	TK
EPHA2 intracellular module	TK
EPHA3	TK

Kinase	Group
EPHA3 intracellular module	TK
EPHA4	TK
EPHA4 intracellular module	TK
EPHA5	TK
EPHA5 intracellular module	TK
EPHA6	TK
EPHA6 intracellular module	TK
EPHA7	TK
EPHA7 intracellular module	TK
EPHA8	TK
EPHA8 intracellular module	TK
EPHB2	TK
EPHB3	TK
EPHB3 intracellular module	TK
EPHB4	TK
EPHB4 intracellular module	TK
EPHB6	TK
ERBB2	TK
ERBB2 intracellular module	TK
ERBB3	TK
ERBB4 intracellular module	TK
FAK	TK
FAK full-length	TK
FER	TK
FES	TK

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

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

KinaseSeeker™ TK Panel (126 Wild-Type and Full-Length Kinases)

Kinase	Group
FGFR2	TK
FGFR2 intracellular module	TK
FGFR3	TK
FGFR3 intracellular module	TK
FGFR4	TK
FGR	TK
FGR full-length	TK
FLT1	TK
FLT1 intracellular module	TK
FLT2	TK
FLT2 intracellular module	TK
FLT3	TK
FLT4	TK
FLT4 intracellular module	TK
FRK	TK
FRK full-length	TK
FYN	TK
FYN full-length	TK
HCK	TK
HCK full-length isoform1	TK
HCK full-length isoform2	TK
IGF1R	TK
IGF1R intracellular module	TK
INSR	TK
INSR intracellular module	TK

Kinase	Group
INSRR	TK
ITK	TK
JAK2	TK
JAK3	TK
KIT	TK
KIT intracellular module	TK
LCK	TK
LCK full-length	TK
LYN	TK
LYN full-length	TK
MET	TK
MET intracellular module	TK
MST1R/RON	TK
MUSK	TK
MUSK intracellular module	TK
NTRK1	TK
NTRK2	TK
NTRK2 intracellular module	TK
NTRK3	TK
NTRK3 intracellular module	TK
PDGFRA	TK
PDGFRA intracellular module	TK
PDGFRB	TK
PDGFRB intracellular module	TK
PTK6	TK

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

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

KinaseSeeker™ TK Panel (126 Wild-Type and Full-Length Kinases)

Kinase	Group
PTK6 full-length	TK
PYK2	TK
PYK2 full-length	TK
RET	TK
RET intracellular module	TK
ROS1	TK
ROS1 intracellular module	TK
SRC	TK
SRC full-length	TK
SRMS	TK
SYK	TK
TEC	TK
TIE1	TK
TIE1 intracellular module	TK
TIE2	TK
TIE2 intracellular module	TK
TNK1	TK
TNK1 full-length	TK
TNK2	TK
TXK	TK
TYRO3	TK
TYRO3 intracellular module	TK
VEGFR2	TK
VEGFR2 intracellular module	TK
YES1	TK

Kinase	Group
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