

**KinaseSeeker-TE™ Panel**

Kinase	Group
ABL1	TK
ABL2	TK
AMPK1	CAMK
AXL	TK
BTK	TK
CAMK2A	CAMK
CAMK2B	CAMK
CAMK2D	CAMK
CAMKK1	Other
CAMKK2	Other
CDK19	CMGC
EGFR	TK
EPHA3	TK
EPHA4 intracellular module	TK
EPHA5	TK
EPHA7	TK
EPHA8	TK
EPHB4 intracellular module	TK
FAK	TK
FER	TK
FGR	TK
FYN	TK
FYN full length	TK
GRK7 full length	AGC
HCK	TK

Kinase	Group
LYN	TK
MLK1/MAP3K9	TKL
MLK3/MAP3K11	TKL
MST1R/RON	TK
MUSK	TK
PDGFRA intracellular module	TK
PHKG1	CAMK
PKG1	AGC
RET	TK
SNARK/NUAK2	CAMK
SRC full length	TK
VEGFR2	TK

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

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

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