

KinaseSeeker-TE™ Panel

Kinase	Group
ABL1	TK
ABL2	TK
AXL	TK
BTK	TK
CAMK2A	CAMK
CAMK2B	CAMK
CAMK2D	CAMK
CAMKK1	Other
CDK19	CMGC
EGFR	TK
EPHA3	TK
EPHA5	TK
EPHA7	TK
EPHA8	TK
FAK	TK
FER	TK
FGR	TK
FYN	TK
HCK	TK
MLK1/MAP3K9	TKL
MLK3/MAP3K11	TKL
MST1R/RON	TK
MUSK	TK
PDGFRA intracellular module	TK
PHKG1	CAMK

Kinase	Group
RET	TK
VEGFR2	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