

## KinaseSeeker™ Complete Panel (415 Kinases)

Kinase	Group
AAK1	Other
ABL1	TK
ABL1 full-length	TK
ABL1 (E255K)	TK
ABL1 (E255V)	TK
ABL1 (F317I)	TK
ABL1 (F317L)	TK
ABL1 (F359V)	TK
ABL1 (G250E)	TK
ABL1 (H396P)	TK
ABL1 (H396R)	TK
ABL1 (L248R)	TK
ABL1 (L248V)	TK
ABL1 (L387M)	TK
ABL1 (M351T)	TK
ABL1 (Q252H)	TK
ABL1 (Q252R)	TK
ABL1 (T315F)	TK
ABL1 (T315I)	TK
ABL1 (T315N)	TK
ABL1 (Y253F)	TK
ABL1 (Y253H)	TK
ABL2	TK
ACVR1	TKL
ACVR1 intracellular module	TKL

Kinase	Group
ACVR1B	TKL
ACVR2A	TKL
ACVR2B	TKL
AKT1	AGC
AKT1 full-length	AGC
AKT2	AGC
AKT2 full-length	AGC
AKT2 (S474A)	AGC
AKT2 (S474D)	AGC
AKT2 (T309A, S474A)	AGC
AKT2 (T309D, S474D)	AGC
AKT3	AGC
AKT3 full-length	AGC
ALK	TK
AMPK	CAMK
AMPK2	CAMK
AMPKa1 full-length	CAMK
AURKA	Other
AURKB	Other
AURKC	Other
AXL	TK
BIKE	Other
BLK	TK
BLK full-length	TK
BMX	TK

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

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

**KinaseSeeker™ Complete Panel (415 Kinases)**

Kinase	Group
BRAF	TKL
BRAF (T599E, S602D)	TKL
BRAF (T599E, V600E, S602D)	TKL
BRAF (V600E)	TKL
BRSK2	CAMK
BTK	TK
BTK full-length	TK
CAMK1	CAMK
CAMK1D	CAMK
CAMK1D full-length	CAMK
CAMK1G	CAMK
CAMK1G full-length	CAMK
CAMK2A	CAMK
CAMK2A full-length	CAMK
CAMK2B	CAMK
CAMK2B full-length	CAMK
CAMK2D	CAMK
CAMK2D full-length	CAMK
CAMK2G	CAMK
CAMK2G full-length	CAMK
CAMKK1	Other
CAMKK1 full-length	Other
CAMKK2	Other
CAMKK2 full-length	Other
CASK	CAMK

Kinase	Group
CASK full-length	CAMK
CDK1	CMGC
CDK19	CMGC
CDK2 - Cyclin E1	CMGC
CDK3	CMGC
CDK4	CMGC
CDK5	CMGC
CDK6	CMGC
CDK7	CMGC
CDK8	CMGC
CDK9	CMGC
CDKL2	CMGC
CDKL3	CMGC
CDKL5	CMGC
CDKL5 full-length	CMGC
CHK1	CAMK
CHK2	CAMK
CK1d	CK1
CK1d full-length	CK1
CK2a2	Other
CLK1	CMGC
CLK1 full-length	CMGC
CLK2	CMGC
CLK2 full-length	CMGC
CLK3	CMGC

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

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

**KinaseSeeker™ Complete Panel (415 Kinases)**

Kinase	Group
CLK3 full-length	CMGC
CSF1R	TK
CSF1R intracellular module	TK
CSK	TK
CSK full-length	TK
DAPK1	CAMK
DAPK1 full-length	CAMK
DAPK2	CAMK
DAPK3	CAMK
DCLK2	CAMK
DDR1	TK
DDR1 intracellular module	TK
DDR2	TK
DDR2 intracellular module	TK
DMPK	AGC
DRAK1	CAMK
DRAK2	CAMK
DYRK1A	CMGC
EGFR	TK
EGFR intracellular module	TK
EGFR (G719C)	TK
EGFR (G719S)	TK
EGFR (L858R)	TK
EGFR (L858R) intracellular module	TK
EGFR (L861Q)	TK

Kinase	Group
EGFR del L747-P753 Ins S	TK
EPHA1	TK
EPHA1 intracellular module	TK
EPHA2	TK
EPHA2 intracellular module	TK
EPHA3	TK
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

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

## KinaseSeeker™ Complete Panel (415 Kinases)

Kinase	Group
ERBB3	TK
ERBB4 intracellular module	TK
ERK5	CMGC
FAK	TK
FAK full-length	TK
FER	TK
FES	TK
FGFR2	TK
FGFR2 intracellular module	TK
FGFR3	TK
FGFR3 intracellular module	TK
FGFR3 (G697C)	TK
FGFR4	TK
FGR	TK
FGR full-length	TK
FLT1	TK
FLT1 intracellular module	TK
FLT2	TK
FLT2 intracellular module	TK
FLT3	TK
FLT3 (D835H)	TK
FLT3 (D835Y)	TK
FLT3 (ITD)	TK
FLT3 (K663Q)	TK
FLT3 (N841I)	TK

Kinase	Group
FLT3 (R834Q)	TK
FLT4	TK
FLT4 intracellular module	TK
FRK	TK
FRK full-length	TK
FYN	TK
FYN full-length	TK
GCN2	STE
GRK7	AGC
GRK7 full-length	AGC
GSK3A	CMGC
GSK3A full-length	CMGC
GSK3B	CMGC
HCK	TK
HCK full-length isoform1	TK
HCK full-length isoform2	TK
HIPK4	CMGC
HIPK4 full-length	CMGC
IGF1R	TK
IGF1R intracellular module	TK
IKKe	Other
IKKe full-length	Other
INSR	TK
INSR intracellular module	TK
INSRR	TK

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

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

## KinaseSeeker™ Complete Panel (415 Kinases)

Kinase	Group
IRAK3	TKL
IRAK3 full-length	TKL
ITK	TK
JAK2	TK
JAK3	TK
KIT	TK
KIT intracellular module	TK
KIT (A829P) intracellular module	TK
KIT (D816H) intracellular module	TK
KIT (D816V) intracellular module	TK
KIT (D820E) intracellular module	TK
KIT (D820Y) intracellular module	TK
KIT (L576P) intracellular module	TK
KIT (N822K) intracellular module	TK
KIT (T670E) intracellular module	TK
KIT (T670I) intracellular module	TK
KIT (V559D) intracellular module	TK
KIT (V654A) intracellular module	TK
KIT (Y823D) intracellular module	TK
LCK	TK
LCK full-length	TK
LIMK1	TKL
LYN	TK
LYN full-length	TK
MAK	CMGC

Kinase	Group
MAP3K10	TKL
MAP3K11	TKL
MAP3K2	STE
MAP3K9	TKL
MAP4K1	STE
MAP4K2	STE
MAP4K2 full-length	STE
MAP4K3	STE
MAP4K5	STE
MAP4K5 full-length	STE
MAP4K6	STE
MAPK11	CMGC
MAPK12	CMGC
MAPK13	CMGC
MAPK14	CMGC
MAPK9 full-length	CMGC
MARK1	CAMK
MARK1 full-length	CAMK
MARK2	CAMK
MARK2 full-length	CAMK
MARK3	CAMK
MARK4	CAMK
MELK	CAMK
MET	TK
MET intracellular module	TK

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

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

## KinaseSeeker™ Complete Panel (415 Kinases)

Kinase	Group
MET (M1250T)	TK
MET (Y1235D)	TK
MKNK2	CAMK
MSK1	AGC
MSK1 full-length	AGC
MSK2	AGC
MSK2 full-length	AGC
MST1	STE
MST1 full-length	STE
MST1R/RON	TK
MST3	STE
MST3 full-length	STE
MUSK	TK
MUSK intracellular module	TK
MYLK	CAMK
MYLK2	CAMK
MYLK2 full-length	CAMK
MYLK4	CAMK
MYLK4 full-length	CAMK
MYT1	Other
MYT1 full-length	Other
NEK5	Other
NEK9	Other
NTRK1	TK
NTRK2	TK

Kinase	Group
NTRK2 intracellular module	TK
NTRK3	TK
NTRK3 intracellular module	TK
NUAK1	CAMK
NUAK1 full-length	CAMK
PAK4	STE
PAK7	STE
PCTK1	CMGC
PCTK2	CMGC
PDGFRA	TK
PDGFRA intracellular module	TK
PDGFRA (D842V)	TK
PDGFRA (D846Y)	TK
PDGFRA (T674I)	TK
PDGFRB	TK
PDGFRB intracellular module	TK
PDK1	AGC
PDK1 full-length	AGC
PfARK1	<i>P. falciparum</i>
PfCDPK2	<i>P. falciparum</i>
PfCDPK5	<i>P. falciparum</i>
PfGSK3	<i>P. falciparum</i>
PfNEK1	<i>P. falciparum</i>
PfNEK3	<i>P. falciparum</i>
PfPK5	<i>P. falciparum</i>

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

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

## KinaseSeeker™ Complete Panel (415 Kinases)

Kinase	Group
PfPK6	<i>P. falciparum</i>
PfPKB	<i>P. falciparum</i>
PFTAIRE1	CMGC
PFTAIRE1 full-length	CMGC
PFTAIRE2	CMGC
PHKG1	CAMK
PIK3CA intracellular module	Lipid
PIK3CA (C420R) intracellular module	Lipid
PIK3CA (E542K) intracellular module	Lipid
PIK3CA (E545A) intracellular module	Lipid
PIK3CA (E545K) intracellular module	Lipid
PIK3CA (Q546K) intracellular module	Lipid
PIK3CA (I800L) intracellular module	Lipid
PIK3CA (M1043I) intracellular module	Lipid
PIK3CA (H1047L) intracellular module	Lipid
PIK3CB intracellular module	Lipid
PIK3CD intracellular module	Lipid
PIK3CG	Lipid
PIM1	CAMK
PIM2	CAMK
PKA	AGC
PKAb	AGC
PKCd	AGC
PKCe	AGC
PKCg	AGC

Kinase	Group
PKCh	AGC
PKCt	AGC
PKD1	CAMK
PKD2	CAMK
PKD3	CAMK
PKG1	AGC
PKG1 full-length	AGC
PKG2	AGC
PKN3	AGC
PKX1	AGC
PLK1	Other
PLK2	Other
PLK3	Other
PLK4	Other
PLK4 full-length	Other
PTK6	TK
PTK6 full-length	TK
PYK2	TK
PYK2 full-length	TK
RET	TK
RET intracellular module	TK
RET (M918T)	TK
RET (V804L)	TK
RET (V804M)	TK
RIPK2	TKL

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

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

## KinaseSeeker™ Complete Panel (415 Kinases)

Kinase	Group
RIPK2 full-length	TKL
ROS1	TK
ROS1 intracellular module	TK
RSK1	AGC
RSK1 full-length	AGC
RSK2	AGC
RSK2 full-length	AGC
RSK3	AGC
RSK3 full-length	AGC
RSK4	AGC
RSK4 full-length	AGC
SGK2	AGC
SGK3	AGC
SIK1	CAMK
SIK1 full-length	CAMK
SIK2	CAMK
SIK2 full-length	CAMK
SIK3	CAMK
SLK	STE
SLK full-length	STE
SNARK	CAMK
SNARK full-length	CAMK
SRC	TK
SRC full-length	TK
SRC (Y530F) full-length	TK

Kinase	Group
SRMS	TK
STK10	STE
STK10 full-length	STE
STK16	Other
STK33	CAMK
STK33 full-length	CAMK
STK35	Other
STK36	Other
SYK	TK
TAOK1	STE
TAOK2	STE
TAOK2 full-length	STE
TAOK3	STE
TBK1	Other
TBK1 full-length	Other
TEC	TK
TESK1	TKL
TESK2	TKL
TESK2 full-length	TKL
TIE1	TK
TIE1 intracellular module	TK
TIE2	TK
TIE2 intracellular module	TK
TNK1	TK
TNK1 full-length	TK

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

Intracellular Module: Construct contains Cytoplasmic Region in Receptor Tyrosine Kinases

**KinaseSeeker™ Complete Panel (415 Kinases)**

Kinase	Group
TNK2	TK
TNNI3K	TKL
TTK	STE
TXK	TK
TYRO3	TK
TYRO3 intracellular module	TK
ULK2	Other
ULK2 full-length	Other
ULK3	Other
VEGFR2	TK
VEGFR2 intracellular module	TK
YANK2	AGC
YES1	TK
YES1 full-length	TK
YSK1	STE

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