

KinaseSeeker™

Kinase Profiling Service

Luceome Biotechnologies
3938 E. Grant Road, # 229
Tucson, Arizona 85712
Ph: (520) 495-0161
www.luceome.com

Luceome Biotechnologies provides a split luciferase based high throughput assay for profiling of small molecules against a panel of kinases, enabling discovery of effective and safer pharmaceuticals. Our *KinaseSeeker* Profiling Service combines the power of split-protein technology with luminescent detection to allow comprehensive screening of compound libraries at cellular ATP concentrations.

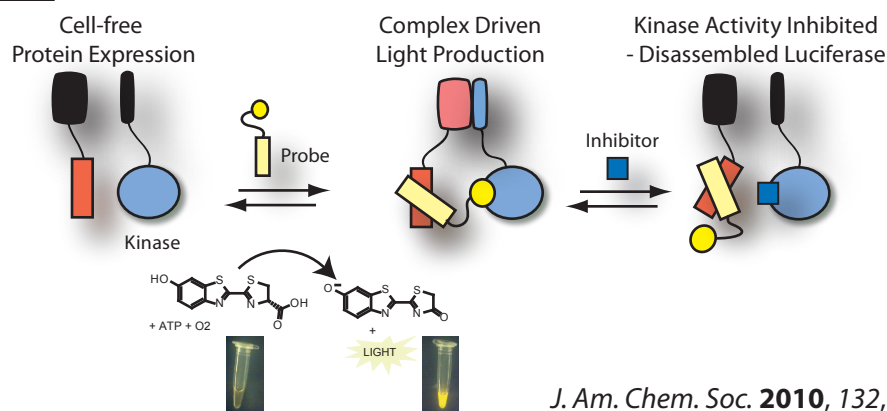
KinaseSeeker™ Technology

KinaseSeeker is a homogeneous binding assay where the displacement of an active site dependent probe by the drug candidate is measured by a change in luminescence signal. Luminescence readout translates into a highly sensitive and robust assay with low background and minimal interference from test compounds.

Features & Benefits

- Luminescence detection: Low false positives from fluorescent compounds
- Binding Assay: Identify active site directed and allosteric inhibitors
- Homogeneous: Portable and adaptable to high-throughput
- Sensitive: Low Background; Broad dynamic range
- Economical: 50 kinase panel profiling - \$400
- Choose the kinases you want
- Fast: 7-10 day turnaround time
- Data kept confidential

ASSAY PRINCIPLE:



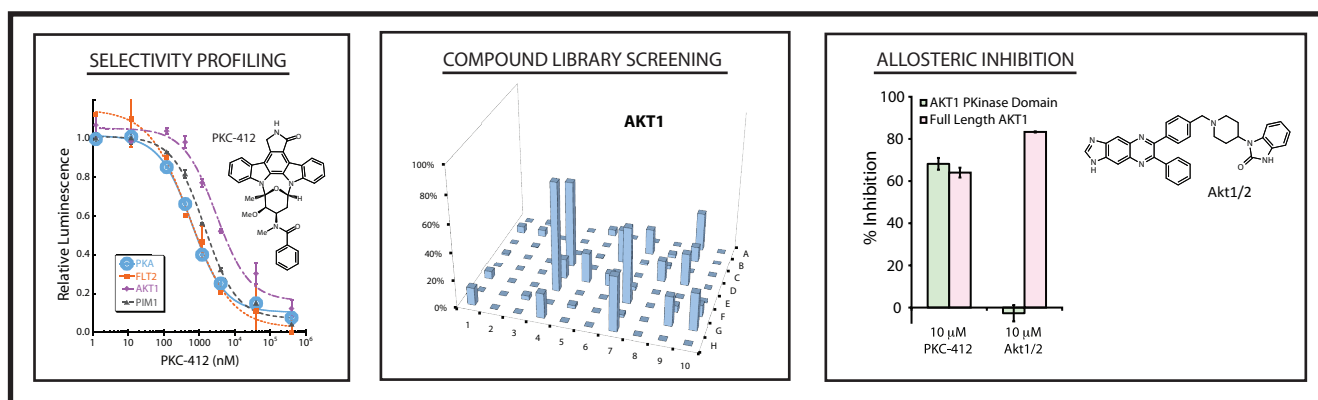
KinaseSeeker™

Kinase Profiling Service

Luceome Biotechnologies
3938 E. Grant Road, # 229
Tucson, Arizona 85712
Ph: (520) 495-0161
www.luceome.com

Available Screening Options:

- (i) Inhibition at single inhibitor concentrations
- (ii) Dose response titrations - (IC_{50}) determination
- (iii) Compound library screening
- (iv) Custom assay development



Available Panels:

- (i) Panel 1: TK-rich (50 kinases, primarily TK/TKL family) - \$400
- (ii) Panel 2: AGC-rich (50 kinases, primarily AGC family) - \$400
- (iii) Panel 3: Mixed panel (50 kinases, representing all families) - \$400
- (iv) Panel 4: Complete panel of 99 kinases - \$750
- (v) A la Carte: \$10/kinase/compound (Minimum order of \$400 required)

Contact:

Reena Zutshi, Ph.D.
Ph: (520) 495-0161
Reena.Zutshi@luceome.com

KinaseSeeker™

Kinase Profiling Service

Luceome Biotechnologies
 3938 E. Grant Road, # 229
 Tucson, Arizona 85712
 Ph: (520) 495-0161
 www.luceome.com

KinaseSeeker™ Panel

ABL2	CAMK2B	IGFR1	PIM1	RPS6KA4
AKT1	CAMK2D	IKK-ε	PIM2	RPS6KA5
AKT1(FL)	CAMKK1	INSR	PKAC-α	RPS6KA6
AKT2	CAMKK2	ITK	PKAC-β	SGK3
AKT2(S474A)	CHEK1	LYN	PKC-δ	SNF1LK
AKT2(S474D)	CLK1	MARK1	PKC-ε	SNF1LK2
AKT2(T309A,S474A)	CLK2	MARK2	PKC-γ	SLK
AKT2(T309D,S474D)	DAPK1	MARK3	PKC-η	SNARK
AKT3	DAPK2	MET	PKC-τ	SRC
AMPK-α1	DDR1	MLK1	PKX	SYK
AMPK-α2	DDR2	MLK3	PRKD2	TBK1
AURKA	DMPK	MST2	PRKD3	TIE1
AURKB	EPHA1	MUSK	PKG1	TNK2
AURKC	FGFR2	p38-γ	PKN3	TRKB
AXL	FLT1	PAK1	PLK4	TRKC
BIKE	FLT2	PAK1(T423A)	PTK2	VEGFR2
BLK	FLT3	PAK1(T423E)	PTK2B	YANK2
CAMK1	FYN	PDGFRA	RPS6KA1	YES1
CAMK1D	GSK3-α	PDGFRB	RPS6KA2	YSK1
CAMK1G	HCK	PDK1	RPS6KA3	