

# KinaseSeeker™

## Kinase Profiling Service

Luceome Biotechnologies  
1665 E. 18th St, Suite 106  
Tucson, Arizona 85719  
Ph: (520) 495-0161  
www.luceome.com

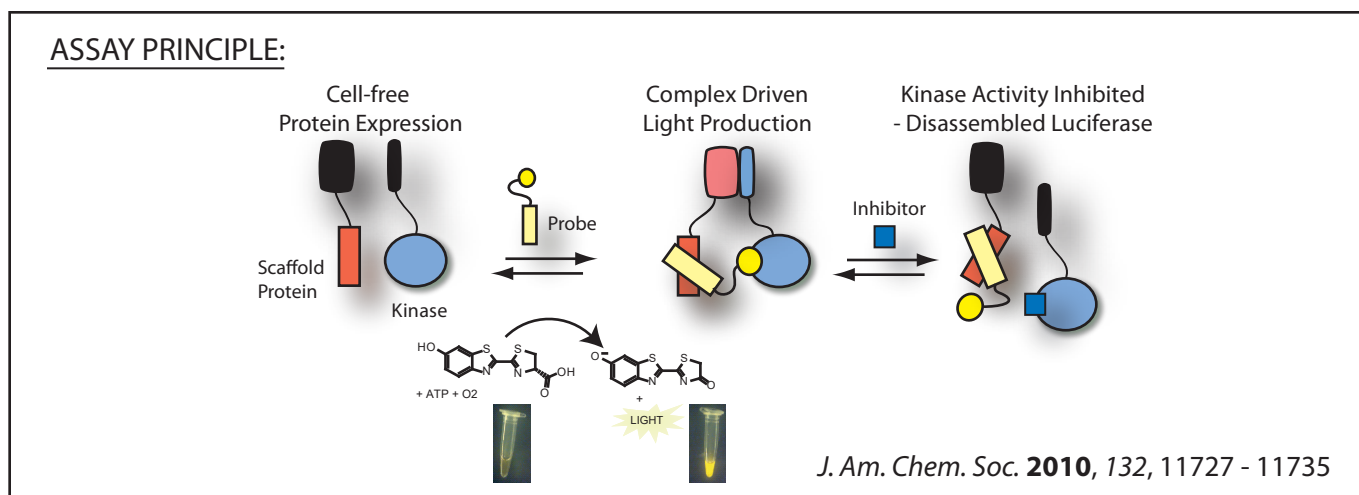
Luceome Biotechnologies provides a split luciferase based high throughput assay for profiling of small molecules against a panel of kinases at cellular ATP concentrations, enabling discovery of effective and safer pharmaceuticals. Our *KinaseSeeker* Profiling Service now includes full-length kinases, enabling discovery of conformation and activation state sensitive kinase inhibitors. The use of full-length kinase constructs also allows identification of allosteric inhibitors that bind the PIF-pocket (PDK1), PH domain (AKTs), SH domains (SFKs), etc.

### **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**

- 415 kinases - Choose from preset kinase panels or build your own panel
- Full-Length Kinases: Evaluate impact of kinase conformation and activation state on compound inhibition
- Binding Assay: Identify active site directed and allosteric inhibitors
- Luminescence detection: Low false positives from fluorescent compounds
- Fast: 10-15 day turnaround time
- Economical: Contact us for pricing ([info@luceome.com](mailto:info@luceome.com))



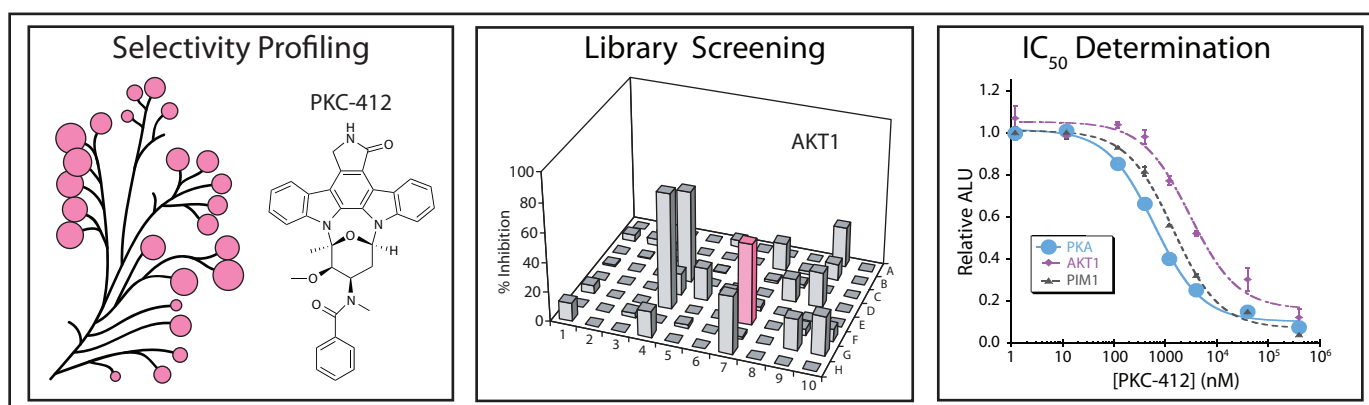
# KinaseSeeker™

## Kinase Profiling Service

Luceome Biotechnologies  
1665 E. 18th St, Suite 106  
Tucson, Arizona 85719  
Ph: (520) 495-0161  
www.luceome.com

### Available Screening Options:

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



Ghosh *et al.*, *J. Am. Chem. Soc.* **2010**, *132*, 11727-11735

Ghosh *et al.*, *J. Med. Chem.* **2012**, *55*, 1526-1537

### Available Panels:

- (i) Full-Length Kinase Panel (112 wild-type full-length kinases distributed across kinome)
- (ii) Tyrosine Kinase Panel (126 wild-type kinase-domain and full-length tyrosine kinases)
- (iii) Mixed panel (50 select kinases distributed across kinome)
- (iv) Plasmodium falciparum Panel (11 *Plasmodium falciparum* kinases)
- (v) Complete panel of 415 kinases
- (vi) A la Carte: Pick kinases of your choice and build your own panel

### Contact:

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