

Novel Gyrasol Sensor Assay Platform with the Hybrid Technology™ Multi-Mode Reader Synergy™ 4

Frauke Rininsland¹, Xavier Amouretti² and Paul Held²

1. Gyrasol Technologies, Santa Fe, NM, 2. BioTek Instruments, Inc., Winooski, VT



Sphingosine Kinase

PI-3 Kinase

PKA - New Platform

Abstract

The Gyrasol Sensor is a novel platform for biochemical kinase / phosphatase assays suitable for small molecule drug discovery. This homogeneous, antibody-free method is universal in nature and is based on fluorescence quench, characterized by signal increase with kinase inhibition. The assay is tolerant to wide concentrations of ATP and substrate, which makes the assay tunable for the inhibition of kinase through either substrate or ATP binding sites and adaptable to either endpoint or kinetic studies. Representative data will be provided for both kinase and phosphatase screening including PI3-K and Sphingosine, noted for their difficult lipid-based substrates. Robust assays, with z' -factors in excess of 0.8 were obtained with the Synergy™ 4 Multi-Mode Microplate Reader using a number of different fluorophores and monochromator settings.

Novel Gyrasol Sensor Assay Platform Overview



- Enzyme activity causes quench of substrate fluorescence
- TAMRA commonly used
- In presence of inhibitors, signal increase assay

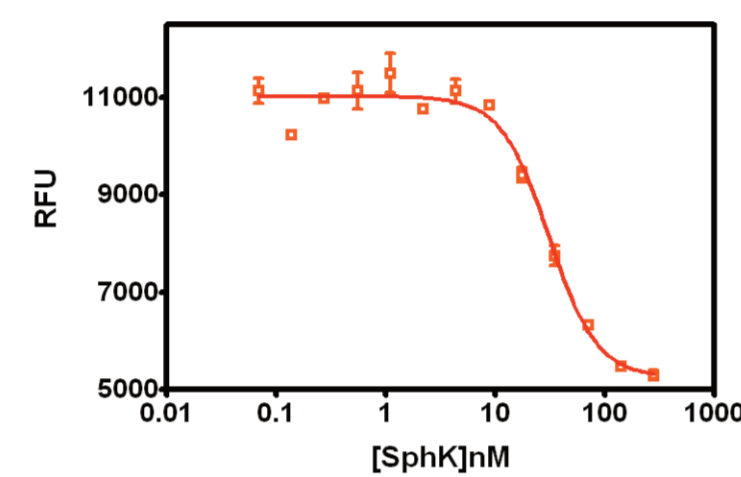
Synergy™ 4 Multi-Mode Reader with Hybrid Technology™



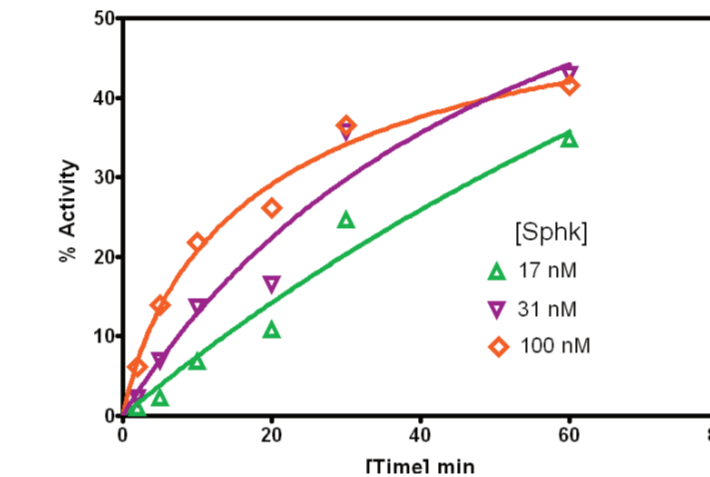
Sphingosine kinase(1) Methods

[Enzyme]: 70 nM (except in titration)
 [Substrate]: 3 μ M
 [ATP]: 100 μ M (+K_m)
 Assay Buffer: Recommended Gyrasol Assay Buffer + 5 mM DTT

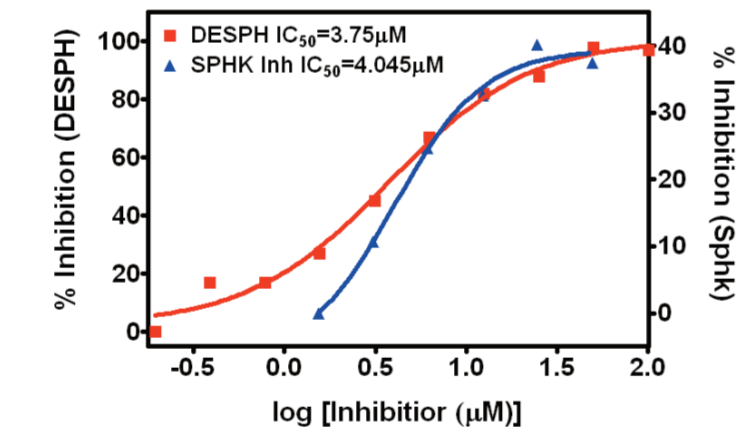
Sphingosine kinase(1) Titration



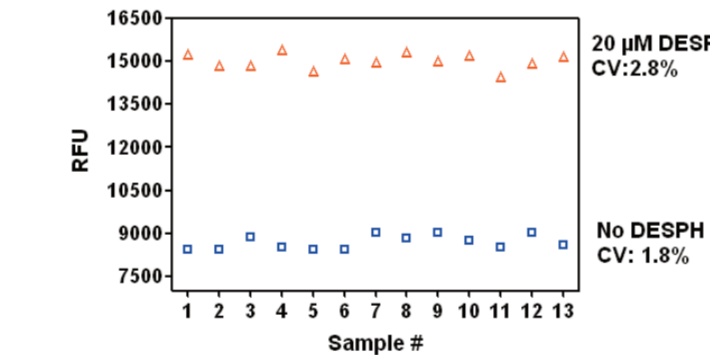
Sphingosine kinase(1) progress curve



Sphingosine kinase(1) Inhibition dose response curves



Sphingosine kinase(1) Statistics



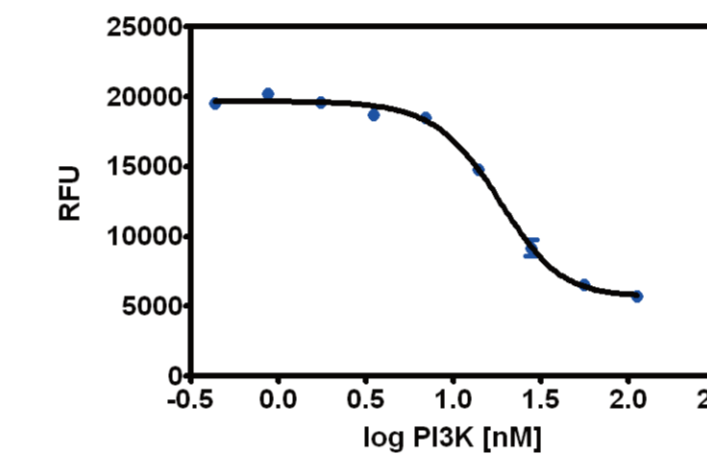
Sphingosine kinase(1) Conclusions

- Sensitive Assays - LOD Sphingosine kinase 7.3 nM
- Determined IC₅₀'s consistent with literature
- Assay Performance suitable for screening applications
 - z' : 0.76
 - CV's < 5%

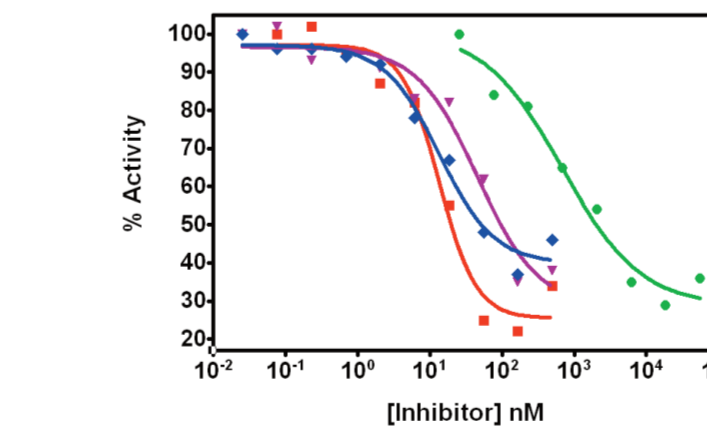
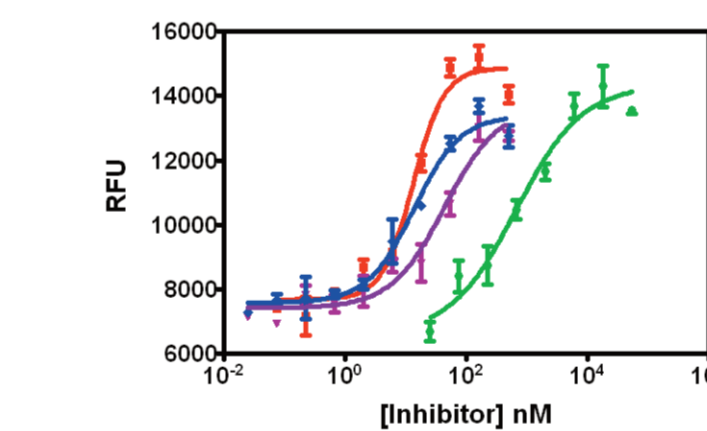
PI-3 Kinase Methods

[Substrate]: 5 μ M
 [ATP]: 150 μ M (+K_m)
 Assay Buffer: Recommended Gyrasol Assay Buffer + 5 mM DTT

PI-3 Kinase Titration



PI-3 Kinase



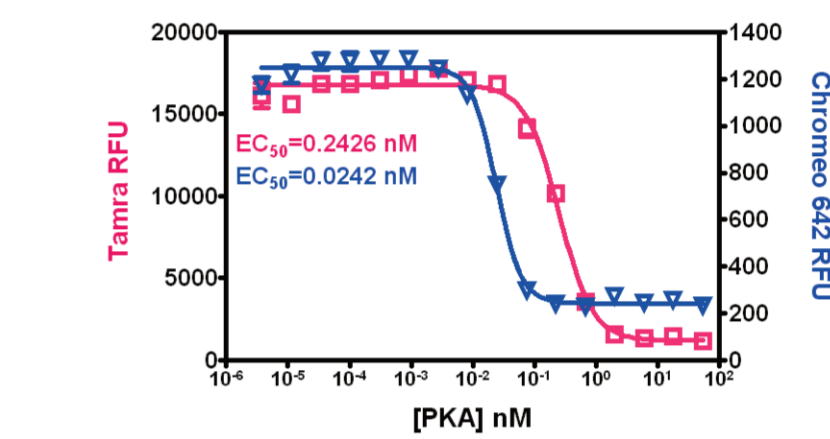
	Gyrasol IC ₅₀ (nM)	Reported IC ₅₀ (nM)
Wortmannin	13	10
PI-103	45	8
PI-3Ka Inhibitor 2	13	2
LY294002	680	1400

PI-3 Kinase Conclusions

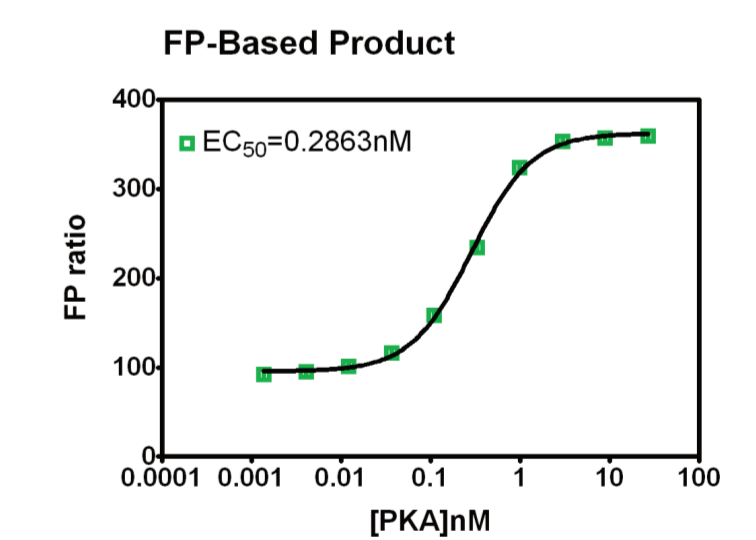
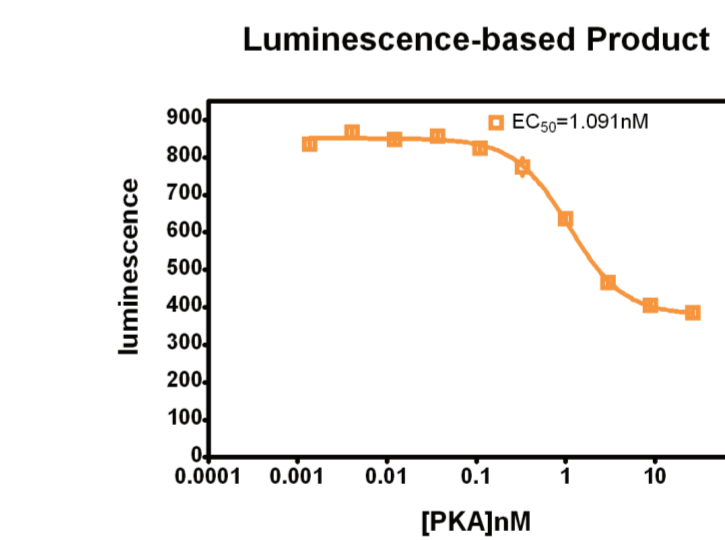
- Sensitive Assays - LOD PI-3 Kinase 1 nM
- Determined IC₅₀'s consistent with literature

New Gyrasol Platform Developments

- Replacement of TAMRA substrate label with Chromeo 642 dye (Active Motif, San Diego)
- Forms complex with the phosphate-associated quencher, thus generating a more favorable distance for electron transfer
- Results in a more sensitive detection



PKA Competitor Comparison with PKA



	LOD (nM)
TAMRA-based Gyrasol	0.2
Chromeo 642-based Gyrasol	0.002
Luminescence-based Product	0.12
FP-based Product	0.005

- Chromeo 642-based Gyrasol is most sensitive