

# Novel Kinase Assays for HTS

Michael Tsionsky,  
Tatiana Plissova, Eli N. Glezer,  
James L. Wilbur, George B. Sigal,  
Hans A. Biebuyck and Jacob N. Wohlstadter



Meso Scale Discovery™  
A division of Meso Scale Diagnostics, LLC.

# Novel Kinase Assays for HTS

## Abstract

Activity assays for Tyrosine Kinase and Serine/Threonine Kinase are highly desirable in drug discovery. There are few satisfactory binding assays for measuring kinase activity, in part because there are few selective and high-affinity antibodies that bind to the phosphorylated products of kinases. Assay methods that use harsh conditions or aggressive wash steps can disrupt weakly bound antibodies and compromise the sensitivity and reliability of the assay. Some homogeneous technologies, which avoid washes, suffer from unacceptable levels of background signal. In this poster, we present kinase assays conducted on a new assay platform developed by Meso Scale Discovery™ (MSD™) that overcomes the difficulties listed above. This platform combines array technologies and electrochemiluminescence detection to achieve sensitive kinase assays that have a simple no-wash format. The assays have been optimized for both 96 and 384 well microplates. Model screens were run on a 10,000 compound library, demonstrating robust assay performance for both Tyrosine and Serine/Threonine Kinases.

### Introduction to Kinases

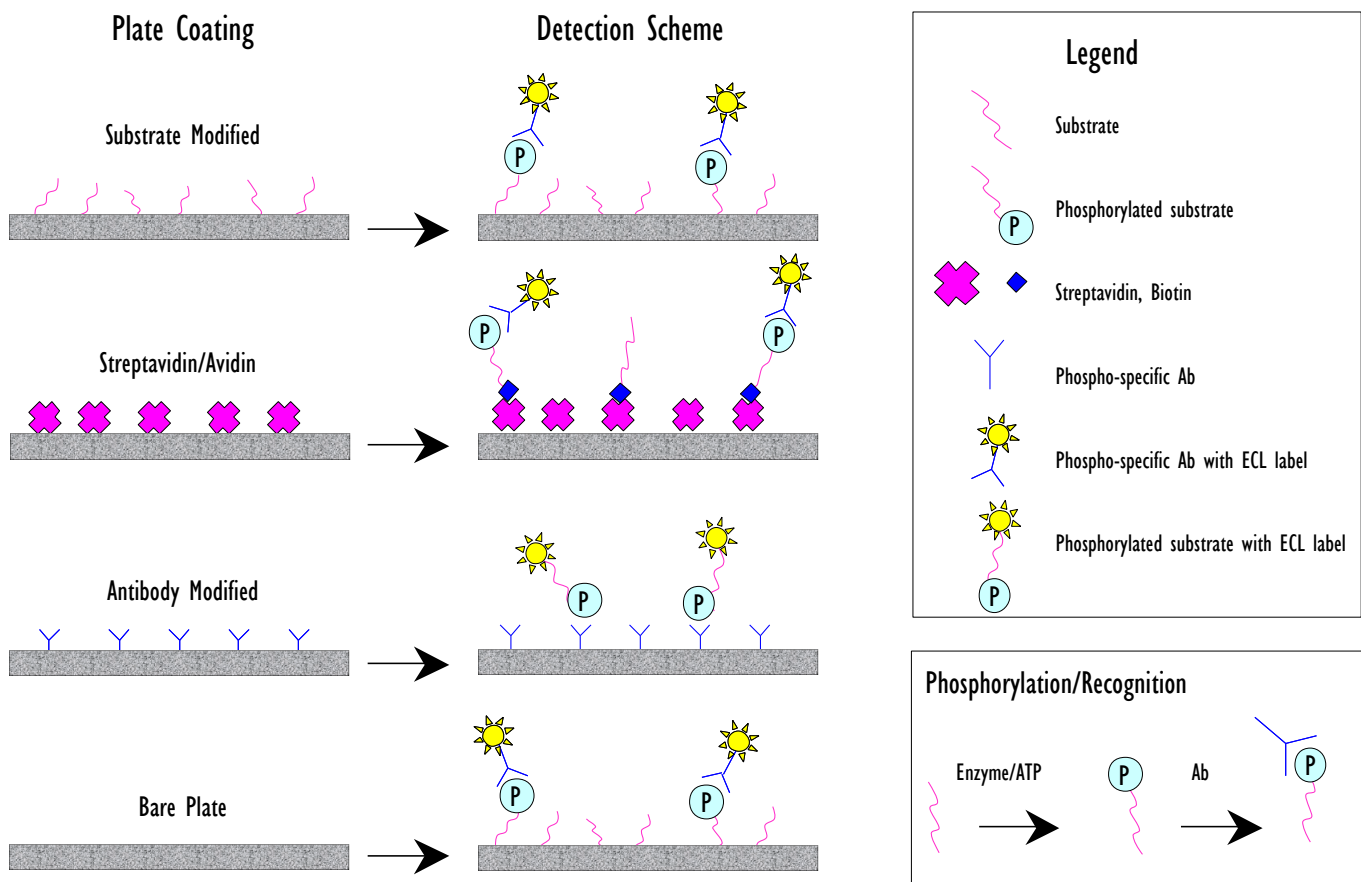
Protein phosphorylation is a critical regulatory mechanism found in cells. Cells often respond to such diverse stimuli as mitogens, inflammatory cytokines, growth factors, and toxins via protein kinases and phosphatases. Modulation of these signaling cascades is often the mechanism for therapies directed against cancer, inflammation, immunology and neurodegenerative disorders<sup>1</sup>. Assays that screen for kinase inhibitors are very popular oncology drug targets<sup>1,2</sup>.



Meso Scale Discovery™  
A division of Meso Scale Diagnostics, LLC.

# Novel Kinase Assays for HTS

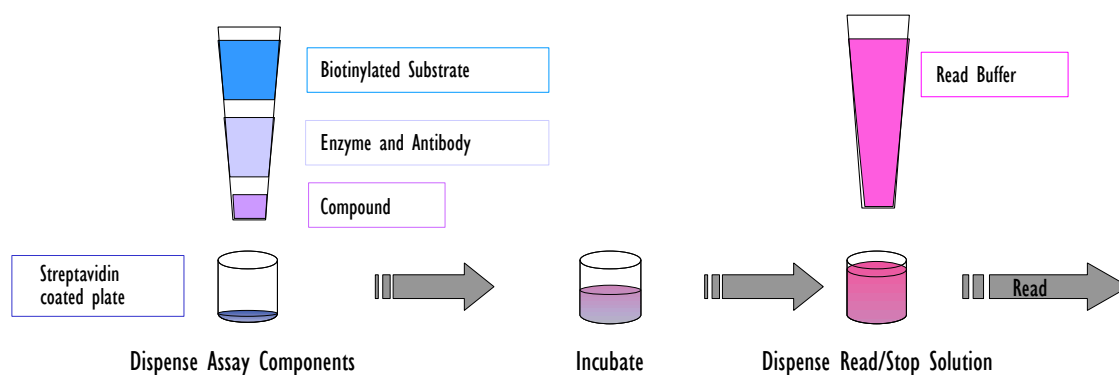
## Versatility of Assay Formats



Meso Scale Discovery<sup>™</sup>  
A division of Meso Scale Diagnostics, LLC.

# Novel Kinase Assays for HTS

## Protocol for Tyrosine Kinase Assay on Streptavidin Plates



### HTS-compatible format

- Only 2 dispense steps
- No wash steps
- Flexible scheduling - only one timed incubation

## Methods

### Plates

Plates were coated by micro-dispensing 2.5  $\mu\text{L}$ /well of 0.5 mg/mL Streptavidin solution using Cartesian's PixSys™ 4200 dispensing system. Plates were then blocked with 5% (w/v) BSA for at least 2 hours. The blocked plates were washed and dried prior to use.

### Assay

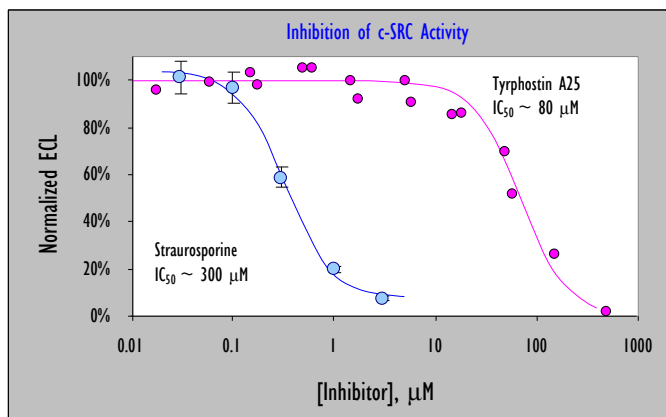
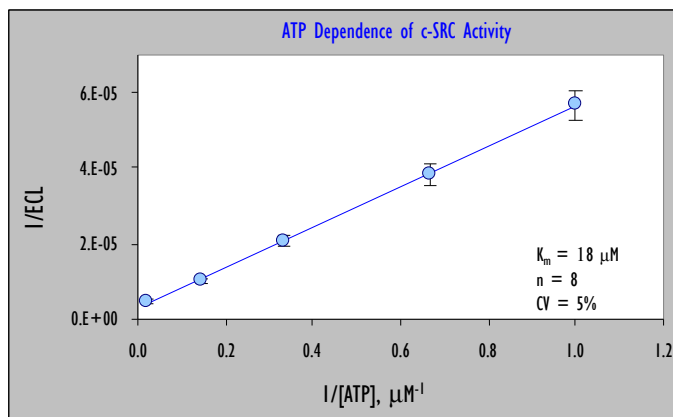
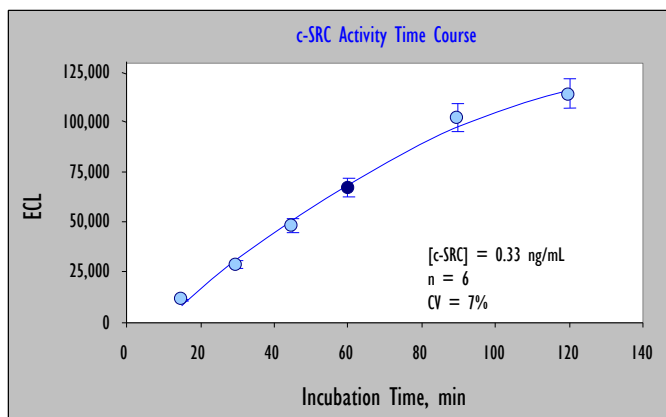
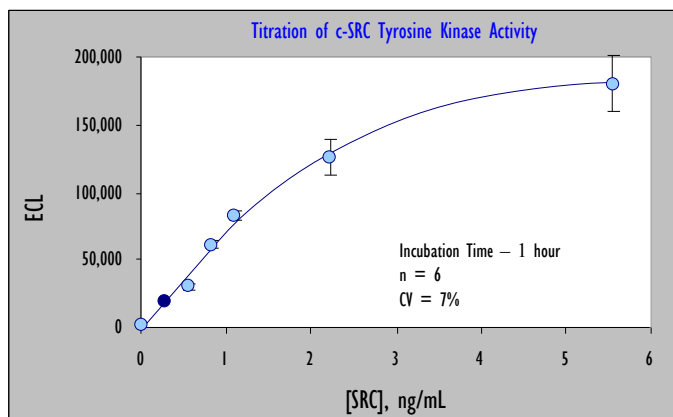
Biotinylated substrate solution (25  $\mu\text{L}$ ), enzyme and labeled Ab solution (25  $\mu\text{L}$ ), and compound from the library (2.5  $\mu\text{L}$ ) were dispensed on Zymark's RapidPlate®. After 1-hour incubation, each well received 200  $\mu\text{L}$  of Read/Stop Buffer solution.



Meso Scale Discovery™  
A division of Meso Scale Diagnostics, LLC.

# Novel Kinase Assays for HTS

## Tyrosine Kinase Assay on 96-Well Substrate Modified Plates



**Selected Conditions:** [c-SRC] = 0.33 ng/mL, [ATP] = 50 mM, labeled Ab = 0.5  $\mu\text{g/mL}$ , poly(Glu,Tyr) 4:1 modified plates, Incubation time = 1 hour

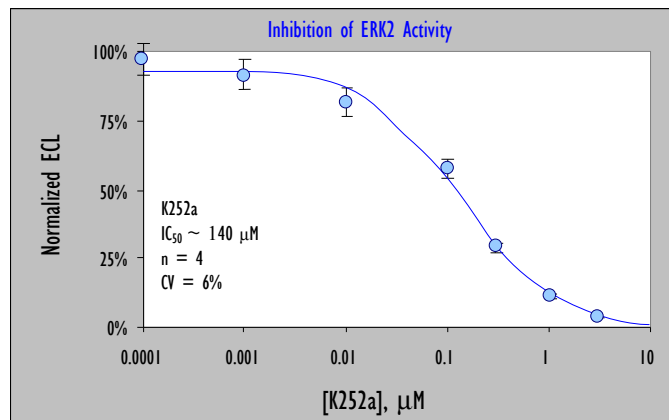
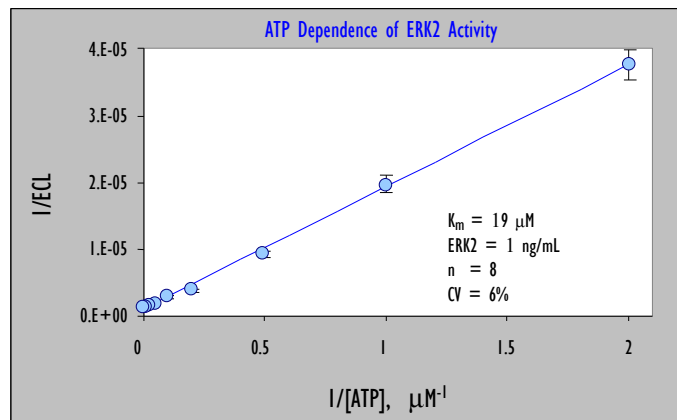
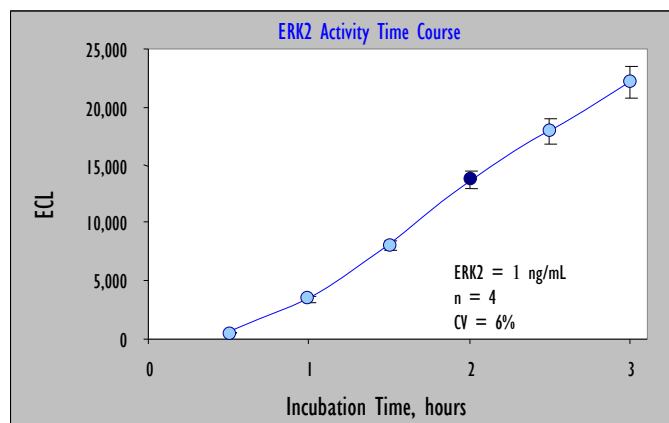
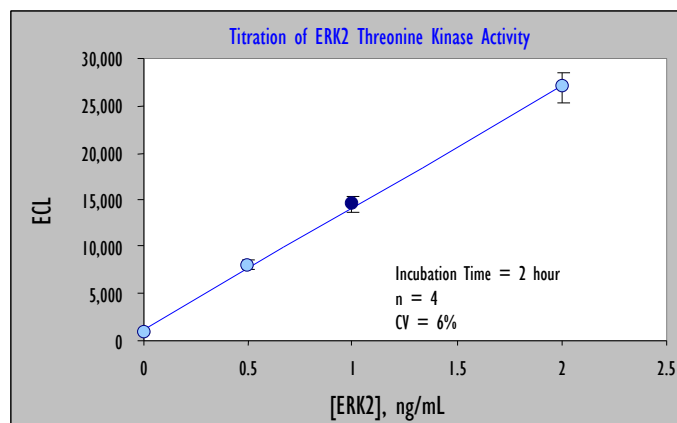
**Assay Performance:** Signal/Background = 60, CV < 10%, Z-factor > 0.6



Meso Scale Discovery<sup>™</sup>  
A division of Meso Scale Diagnostics, LLC.

# Novel Kinase Assays for HTS

## Threonine Kinase Assay on 96-well Streptavidin Modified Plates



**Selected Conditions:** [ERK2] = 1 ng/mL, [ATP] = 50  $\mu\text{M}$ , [bMBP] = 100 nM, primary Ab = 2 nM, secondary Ab = 5 nM, Incubation time = 2 hours

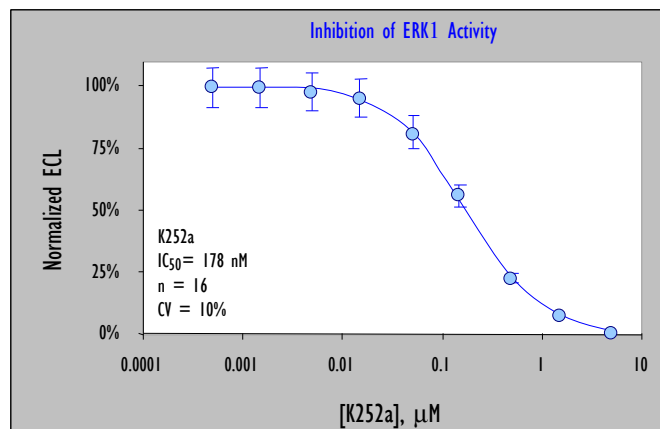
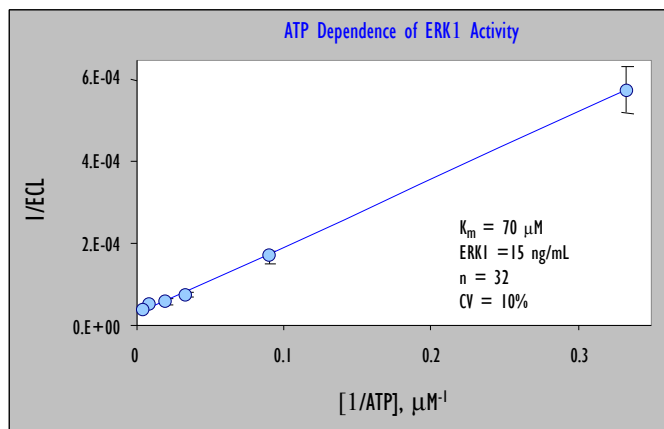
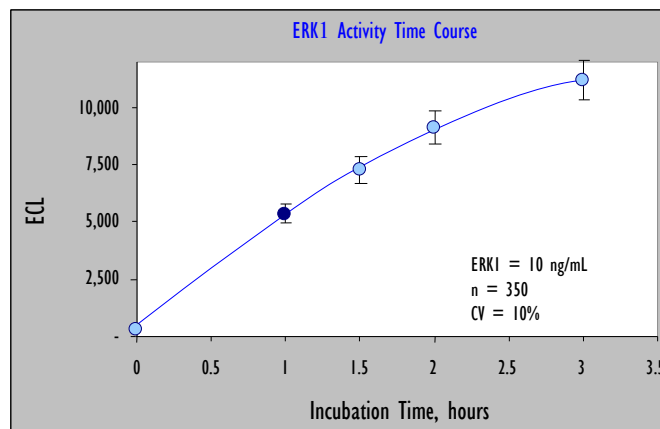
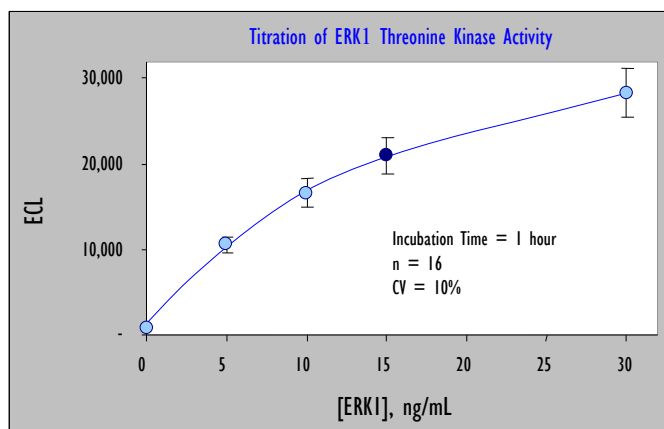
**Assay performance:** Signal/Background = 30, CV < 10%, Z-factor > 0.6



Meso Scale Discovery<sup>™</sup>  
A division of Meso Scale Diagnostics, LLC.

# Novel Kinase Assays for HTS

## Threonine Kinase Assay on 384-well Bare Plates



**Selected Conditions:** [ERK1] = 15 ng/mL, [ATP] = 100  $\mu\text{M}$ , [MBP] = 600 nM, primary Ab = 10 nM, secondary Ab = 25 nM, Incubation Time = 1 hour

**Assay Performance:** Signal/Background = 29, CV = 10%, Z-factor > 0.6

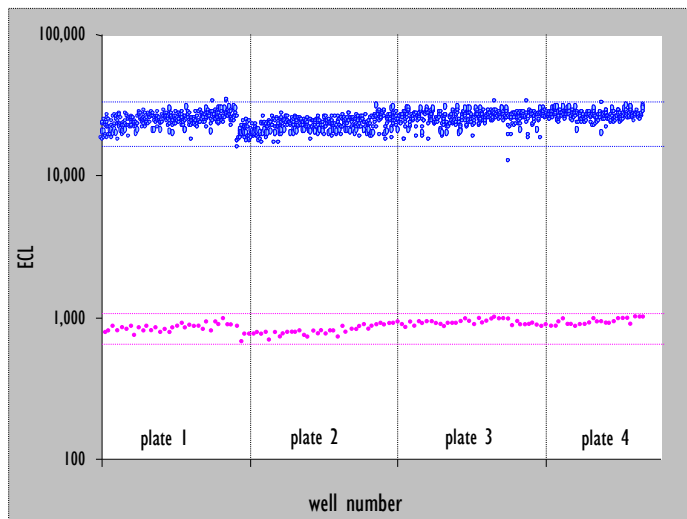


Meso Scale Discovery<sup>™</sup>  
A division of Meso Scale Diagnostics, LLC.

# Novel Kinase Assays for HTS

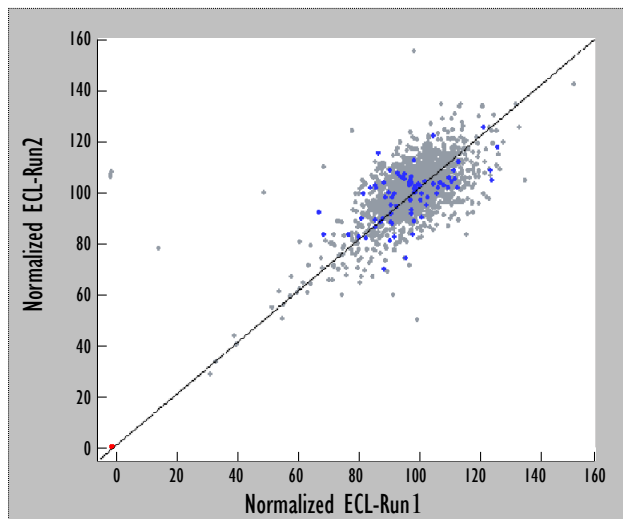
## Assay Performance in HTS applications

Threonine Kinase Controls  
Bare 384-well plates



Signal to Background = 29  
Z-factor = 0.63  
Positive Control CV = 12%  
Negative Control CV = 8%  
Lines represent 3 SD

Tyrosine Kinase Screen in Duplicates  
Substrate-modified 96-well plates



Signal to Background = 60  
Z-factor = 0.75  
Positive Control CV = 8%  
Negative Control CV = 7%  
1,800 compound run shown



Meso Scale Discovery<sup>™</sup>  
A division of Meso Scale Diagnostics, LLC.



# Novel Kinase Assays for HTS

## Conclusion

Tyrosine and Serine/Threonine Kinase assays were developed on a novel platform from Meso Scale Discovery. This platform uses Multi-Array technology that combines array technology with electrochemiluminescent detection.

Multi-Array technology provides a sensitive, robust, simple, non-radioactive means to assay protein kinase activity.

Assay protocols were designed for HTS — no-wash format, short incubation time, automation-friendly protocol.

High quality compound screening data were obtained in 96 and 384-well plates — S/B > 30, CVs — 10-12%, Z-factor > 0.6.

>100-fold reduction in enzyme consumption results from highly sensitive detection (relative to other technologies<sup>3,4</sup>).

Several assay formats were shown to be effective.

### References:

1. Society for Medicines Research Committee (1999). *Drug News Perspect.* 12: 247-251.
2. Levitzki, A. and Gazit, A. (1995). *Science* 267: 1782-1788.
3. Amersham Pharmacia Biotech Catalog, 2000, p.76
4. Packard BioScience-Alpha Screen c-Src Kinase Assay. AN003-ASc Application Note, [www.packardbioscience.com](http://www.packardbioscience.com).



Meso Scale Discovery™  
A division of Meso Scale Diagnostics, LLC.