

# MAP Kinase Activity Assays on MSD Multi-Array™ Platform

Michael Tsionsky  
Tatiana Plissova  
Eli N. Glezer  
and Jacob N. Wohlstadter



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

# MAP Kinase Activity Assays on MSD Multi-Array™ Platform

## 1 Abstract

Mitogen-Activated Protein Kinases (MAPKs) are a widely conserved family of Serine/Threonine protein kinases involved in many cellular programs such as cell proliferation, cell differentiation, cell movement and cell death. MAPK signaling cascades are organized into three-tiered modules. MAPKs are phosphorylated and activated by MAPK-kinases (MAPKK), which in turn are phosphorylated and activated by MAPKK-kinases (MAPKKKs). The MAPKKKs are in turn activated by interaction with a family of small GTPases and/or other protein kinases connecting the MAPK module to the cell surface receptor or external stimulus.

We demonstrate a multiplexed assay approach for monitoring the activity of the entire MAPK cascade or any part of it using the Meso Scale Discovery Multi-Array™ platform. Whole proteins or short synthetic peptides can be employed as substrates. The assay protocols are compatible with high-throughput screening (HTS) and provide sensitive detection of enzyme activity across the MAP Kinase pathways.



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

# MAP Kinase Activity Assays on MSD Multi-Array™ Platform

## 2 Meso Scale Discovery Multi-Array Technology

### Instrument Features

- Highly sensitive
- SECTOR™ Imager 6000 designed for high-throughput screening (HTS)
- SECTOR™ PR 100 Reader ideal for assay development
- Custom optics
- High-speed motion control systems
- Electrochemiluminescence (ECL) detection

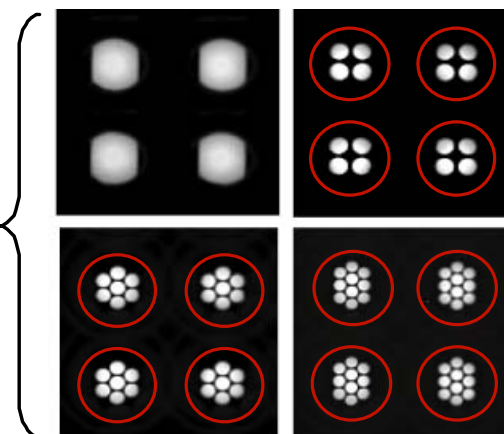
SECTOR™ PR 100 Reader



SECTOR™ Imager 6000

### Plate Features

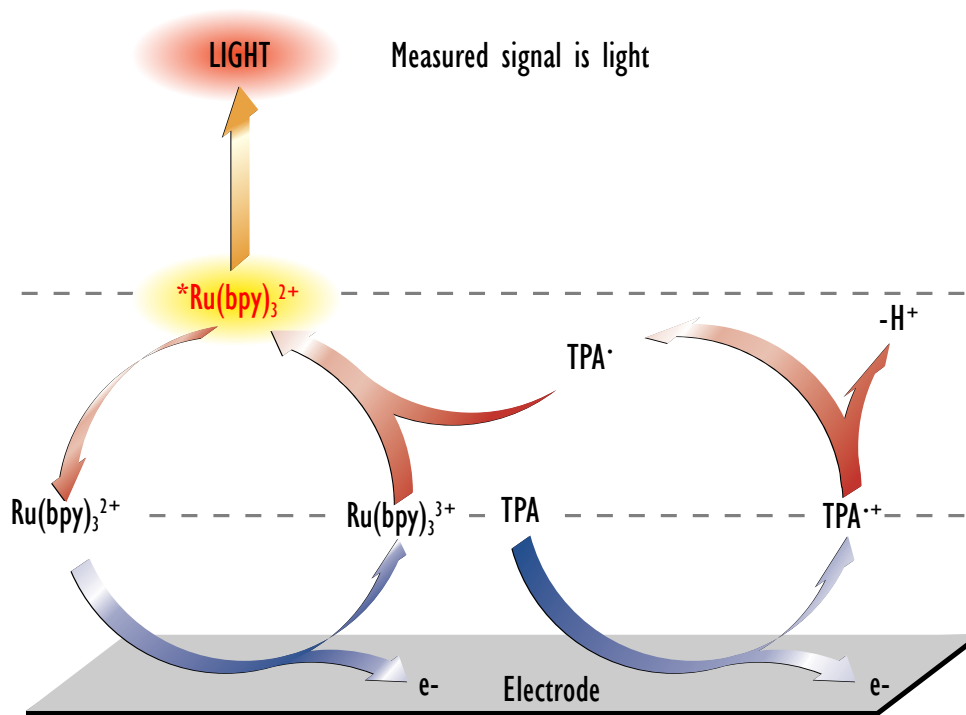
- Disposable Plates
- Carbon Electrodes with high binding capacity
- Suitable electrochemistry for ECL
- Biocompatible: direct immobilization of avidin, IgG, membrane fragments, intact cells, etc.
- Functional Assays: simple binding reactions, GPCRs, enzyme cascades, post-translational modification, etc.



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

# MAP Kinase Activity Assays on MSD Multi-Array™ Platform

## 3 Electrochemiluminescence (ECL)



**Luminescence**

Emitting Light

**Chemi-**

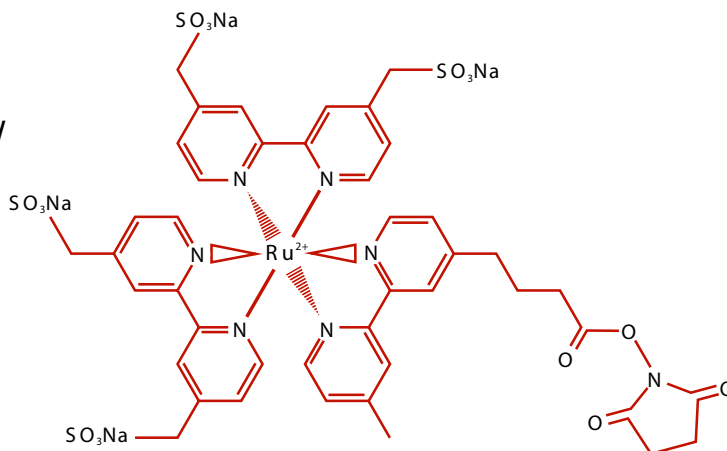
Chemical Energy

**Electro-**

Electrochemically Initiated

### MSD Sulfo-TAG™ Label

- Selective
- Convenient chemistry
- Robust, stable
- Few interferences



- Size, MW: ~1200 daltons
- Stability: Years
- Solubility: Aqueous, DMSO
- Functionality: Hydrophilic
- Specificity: High

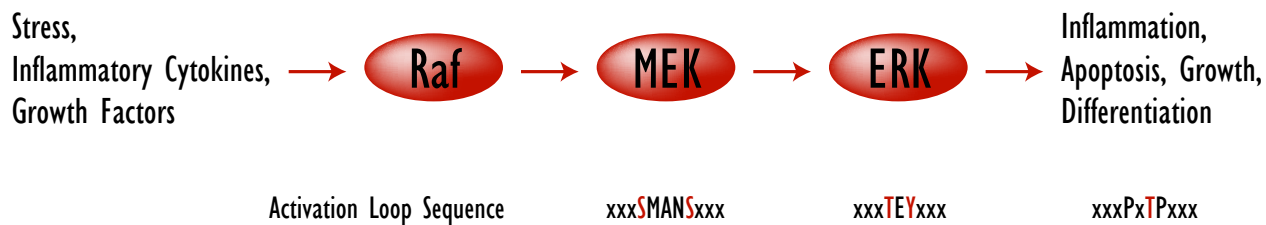


Meso Scale Discovery™  
A division of Meso Scale Diagnostics, LLC.

# MAP Kinase Activity Assays on MSD Multi-Array™ Platform

## 4 MAP Kinase Cascade

### Example of MAPK Cascade



### Substrates Used for MAP Kinase Cascade Signaling

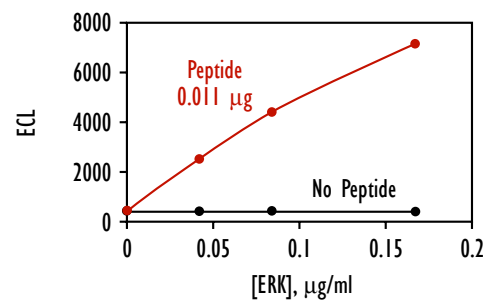
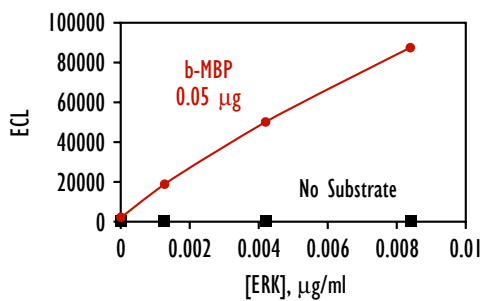
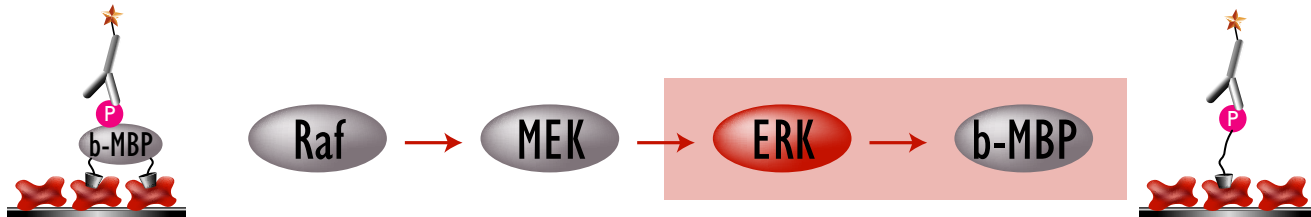
- Whole Proteins: Surface-bound Myelin Basic Protein (MBP), whole ERK and MEK molecules
- Synthetic peptides: Biotinylated short peptides, which contain xxxPxTPxxx, xxxTEYxxx or xxxSMANSxxx motifs, which mimic the target sequences of MBP, ERK and MEK respectively



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

# MAP Kinase Activity Assays on MSD Multi-Array™ Platform

## 5 ERK Activity Assays



**Enzyme:** Active ERK-2

**Substrates:** Biotinylated-MBP or biotinylated synthetic peptide containing P RTP motif immobilized on the surface of MSD Multi-Array 96-well plates coated with streptavidin

**Antibodies:** Mouse anti-phospho-MBP IgG (primary) and MSD Sulfo-TAG™-labeled anti-mouse IgG (secondary)

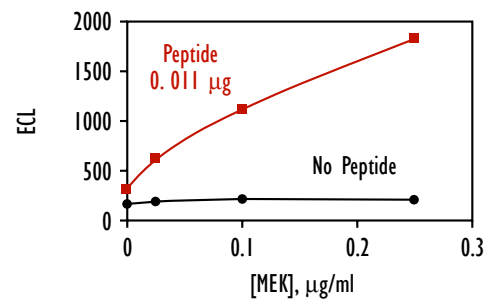
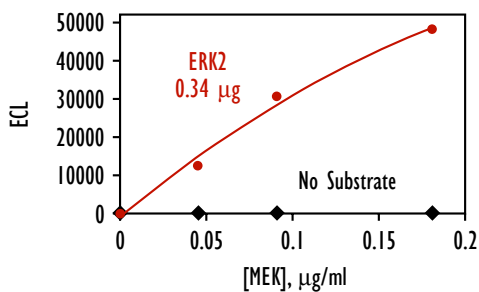
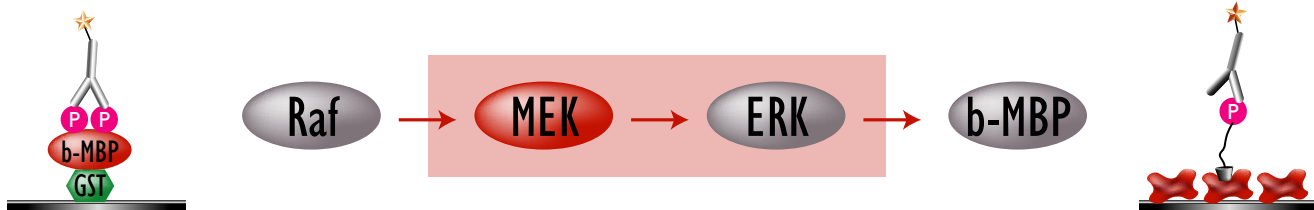
- Assay measures the activity of ERK kinase
- Linear response to ERK concentration
- Low background in the absence of ERK
- Signal-to-background ratio is  $\sim 45$  at 0.001  $\mu\text{g/ml}$  of ERK, with MBP used as the substrate
- Signal-to-background ratio is  $\sim 14$  at 0.1  $\mu\text{g/ml}$  of ERK, with a short peptide substrate



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

# MAP Kinase Activity Assays on MSD Multi-Array™ Platform

## 6 MEK Activity Assays



Enzyme: **Active MEK-1**

Substrates: Inactive GST-ERK-2 or biotinylated synthetic peptide containing **SMANS** motif immobilized on MSD Multi-Array 96-well plates coated with streptavidin

Antibodies: Mp44 Mouse IgG (primary) with MSD Sulfo-TAG-labeled anti-mouse IgG (secondary) or Rp44 Rabbit IgG (primary) with MSD Sulfo-TAG-labeled anti-rabbit IgG (secondary)

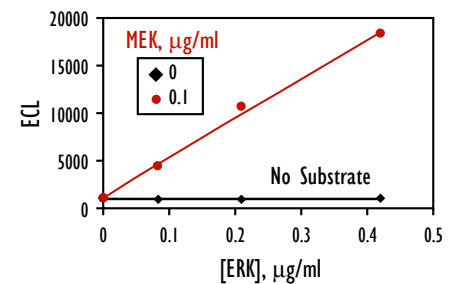
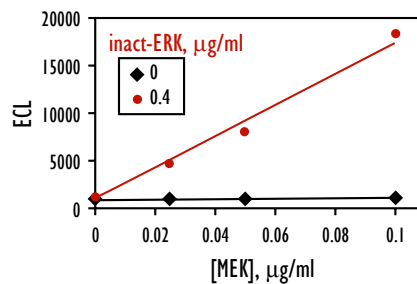
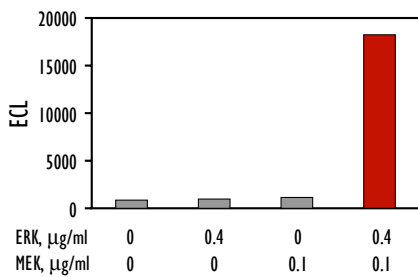
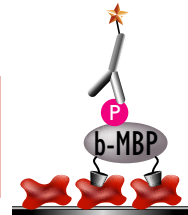
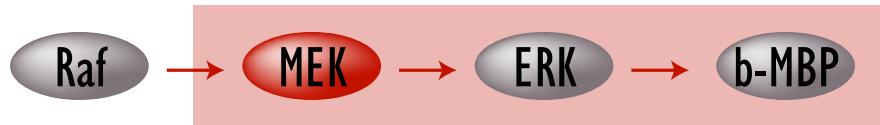
- Assay measures the activity of MEK kinase
- Approx. linear response to MEK concentration
- Low background in absence of MEK
- Signal-to-background ratio at 0.1 µg/ml of MEK is ~80 for whole ERK protein and ~5 for the peptide substrate



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

# MAP Kinase Activity Assays on MSD Multi-Array™ Platform

## 7 Combined MEK & ERK Activity Assay



Enzyme: Active MEK-1, inactive ERK-2

Substrates: Biotinylated-MBP immobilized on MSD Multi-Array  
96-well plates coated with streptavidin

Antibodies: Mouse anti-phospho-MBP IgG (primary),  
MSD Sulfo-TAG-labeled anti-mouse IgG (secondary)

- Assay measures the activity of MEK and ERK kinases simultaneously
- Linear response to both MEK and ERK
- Low background in absence of either enzyme
- Signal-to-background ratio is  $\sim 20$  at 0.4  $\mu\text{g/ml}$  ERK and 0.1  $\mu\text{g/ml}$  MEK

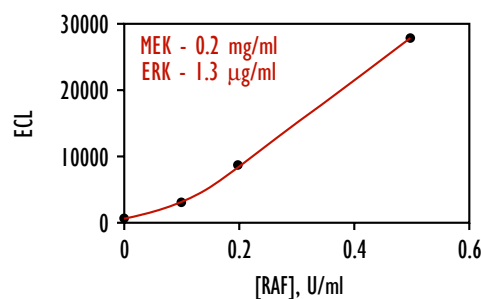
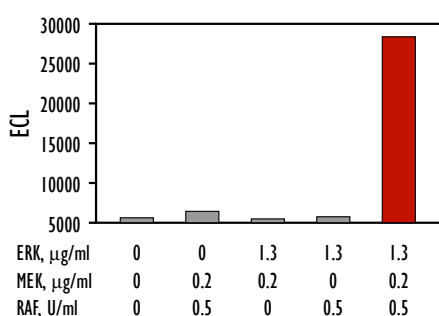
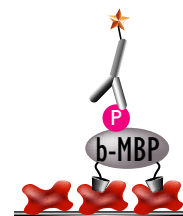
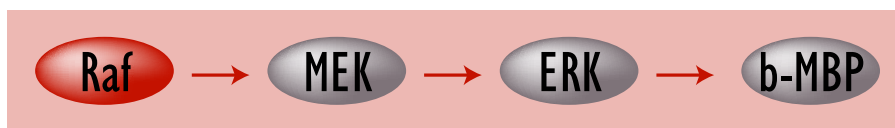


Meso Scale Discovery™  
A division of Meso Scale Diagnostics, LLC.



# MAP Kinase Activity Assays on MSD Multi-Array™ Platform

## 8 Assay for Entire MAPK Cascade



Enzyme: **Active Raf-1**; inactive MEK-1 and ERK-2  
Substrates: Biotinylated-MBP immobilized on MSD Multi-Array  
96-well plates coated with streptavidin  
Antibodies: Mouse anti-phospho-MBP IgG (primary),  
MSD Sulfo-TAG-labeled anti-mouse IgG (secondary)

- Assay simultaneously monitors the activity of entire 3-step cascade
- Linear range exists for all 3 enzymes
- Low background in absence of any components of the MAPK cascade

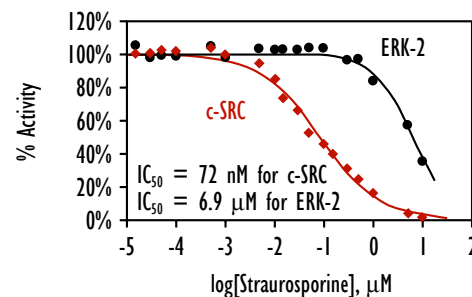
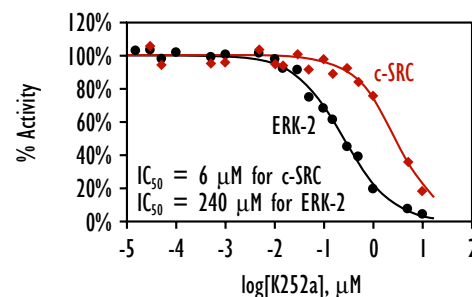
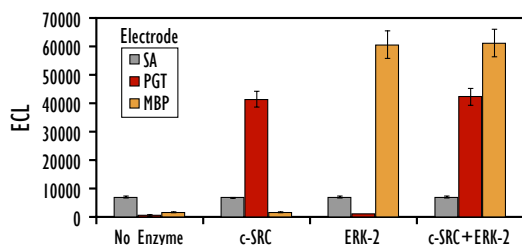
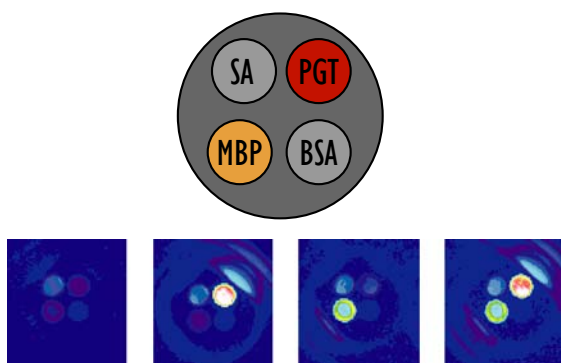


Meso Scale Discovery®  
A division of Meso Scale Diagnostics, LLC.

# MAP Kinase Activity Assays on MSD Multi-Array™ Platform

## 9 Multiplexing: Serine/Threonine and Tyrosine Kinases in one well

Enzyme: Active ERK-2 and Active c-SRC  
Substrates: poly-Glu-Tyr (PGT) and MBP immobilized on the surface of MSD Multi-Spot™ 4-spot plate  
Antibodies: Mouse anti-phospho-MBP IgG, Mouse pY-20, Sulfo-TAG-labeled Anti-Mouse IgG



- Multiplexed format independently monitors the activities of 2 kinases
- Example shows preferential inhibition of ERK-2 by K252a, and c-SRC by Staurosporine

## 10 Conclusions

- Activity of the entire MAPK cascade or any part of it can be monitored using the MSD Multi-Array platform.
- Whole proteins or short synthetic peptides can be employed as substrates.
- Assay protocols are suitable for HTS applications.
- Multiple kinases can be assayed simultaneously in a single sample by using specific target substrates immobilized on a MSD Multi-Spot plate.



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