

## Anti-phospho-MEK1 (Thr292)

Rabbit Polyclonal Antibody

**Catalog # M02-65BR**

Lot # J1274-21

### Cited Applications

WB

Suggested Dilutions:  
WB 1:1,000

Ideal working dilutions for each application should be empirically determined by the investigator.

### Specificity

Recognizes the MEK1 protein phosphorylated at threonine 292

### Cross Reactivity

Human, Mouse, Rat, Bovine, Canine, Chicken and non-Human Primates

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr292 conjugated to KLH

### Formulation

100 µl in 10 mM HEPES (pH 7.5), 150 mM NaCl, 100 µg per ml BSA and 50% glycerol.

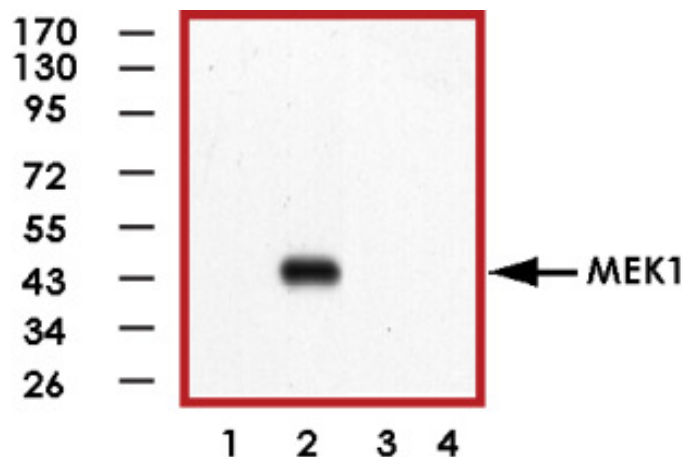
### Scientific Background

MEK1 is a member of the dual specificity protein kinase family and is also referred to as MAP Kinase Kinase or MKK. MEK1 lies upstream of MAPK/ERK and stimulates the enzymatic activity of MAPK/ERK upon a wide variety of extra- and intracellular signals. As an essential component of the MAPK/ERK signal transduction pathway, MEK1 is involved in many cellular processes including proliferation, differentiation, transcriptional regulation and development (1). MEK1 plays a key role in synaptic plasticity in the brain (2). Constitutive activation of MEK1 results in cellular transformation. Thus, MEK 1 represents a likely target for pharmacological intervention in proliferative diseases such as cancer (3). MAP kinase can phosphorylate MEK1 at Thr292 and Thr386.

### References

1. Seger, R. et al: The MAPK signaling cascade. FASEB J. 1995 Jun;9(9):726-35.
2. Adams, J P. et al: Molecular psychology: Roles for the ERK MAP kinase cascade in memory. Annu Rev PharmacolToxicol 2002 42:135-163.
3. Sebolt-Leopold, J S. et al: Blockade of the MAP Kinase Pathway Suppresses growth of colon tumors in vivo. Nature Med. 1999 5: 810-816.

### Sample Data



Western blot of recombinant wild type and mutant MEK1 showing immunolabeling of the ~45kDa MEK1 protein phosphorylated at Thr292. Lanes 1 and 2 are WT MEK1 and Lanes 3 and 4 are mutant MEK1 (T292A). MAP kinase was co-expressed in the samples run in Lanes 2 and 4.

## Anti-phospho-MEK1 (Thr292)

Rabbit Polyclonal Antibody

Catalog Number

M02-65BR

Specific Lot Number

J1274-21

Purification Affinity chromatography

Stability 1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**