

## Anti-MEK1

Mouse Monoclonal Antibody

Catalog # M02-60M

Lot # D316-6

### Cited Applications

- Western blot (1:1000)
- Immunoprecipitation (2 µg)

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

### Specificity

Recognizes the MEK1 protein

### Cross Reactivity

- Western blot human MEK1
- Immunoprecipitation of human MEK1

MEK1 from other species may also be detectable

### Host/Isotype/Clone#

Mouse, IgG1, 7F1

### Immunogen

Recombinant full length human MEK1 protein

### Formulation

0.1M TRIS pH 8.0, 0.1M glycine

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) for several months, and at -20°C for longer periods. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

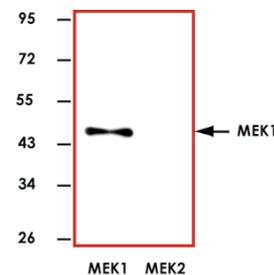
### Scientific Background

MEK1 is a member of the dual specificity protein kinase family that acts as a mitogen-activated protein kinase (MAPK) kinase. 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 MAPK/ERK signal transduction pathway, MEK1 is involved in many cellular processes such as proliferation, differentiation, transcription regulation and development (1). Constitutive activation of MEK1 results in cellular transformation. Thus, MEK1 represents a likely target for pharmacologic intervention in proliferative disease such as cancer (2).

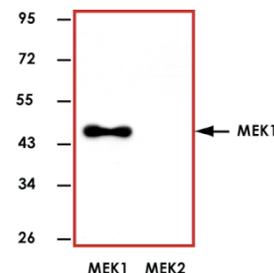
### References

1. Seger, R. et al: The MAPK signaling cascade. FASEB J. 9: 726-735, 1995.
2. Sebolf-Leopold, J.S. et al: Blockade of the MAP kinase pathway suppresses growth of colon tumors in vivo. Nature Med. 5: 810-816, 1999.

### Sample Data



Representative western blot with Anti-MEK1 (1:1000) using 20ng of MEK1 and MEK2 recombinant protein.



Representative immunoprecipitation with Anti-MEK1 (2 µg) using 20ng of MEK1 and MEK2 recombinant protein.

## Anti-MEK1

Mouse Monoclonal Antibody

Catalog Number

M02-60M

Specific Lot Number

D316-6

Purification

Affinity purified using Protein G

Concentration

1.0 µg/µL

Stability

1 yr 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.**