

## MEK1, Unactive

Recombinant full-length human protein expressed in Sf9 cells

**Catalog # M02-14H**

Lot # N053-2

### Product Description

Recombinant human full-length MEK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The MEK1 gene accession number is [NM\\_002755](#).

### Gene Aliases

MAP2K1, MKK1, MAPKK1, PRKMK1

### Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 2mM DTT, 25% glycerol.

### Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

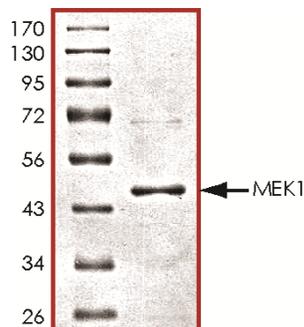
### Scientific Background

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. Sebolt-Leopold, J S. et al: Blockade of the MAP kinase pathway suppresses growth of colon tumors in vivo. *Nature Med.* 5: 810-816, 1999.

### Purity



The purity of unactive MEK1 was determined to be **>90%** by densitometry, approx. MW **47 kDa**.

### Upstream Active Kinases

BRAF, Active	Cat # B08-11G
BRAF(V599E), Active	Cat # B08-13G
COT, Active	Cat # M16-11G
MEK1, Active	Cat # M09-11G
MEK2, Active	Cat # M10-10G
RAF1 (EE), Active	Cat # R01-13G

## MEK1, Unactive

Recombinant full-length human protein expressed in Sf9 cells

Catalog Number M02-14H

Specific Lot Number N053-2

Purity	>90%
Concentration	0.2 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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.**