

## MEK4 (MAP2K4), Unactive

Full-length recombinant human protein expressed in Sf9 cells

Catalog # M05-15G

Lot # K305-1

### Product Description

Recombinant human MEK4 (33-end) was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM\\_003010](#).

### Gene Aliases

MAP2K4, JNKK, MEK4, JNKK1, SERK1, MAPKK4, PRKMK4, SEK1

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 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.

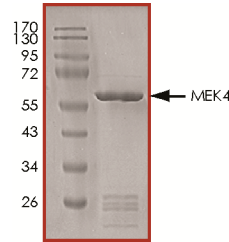
### Scientific Background

MEK4 is a member of the mitogen-activated protein kinase family and is involved in the stress-activated protein kinase pathway where it phosphorylates c-Jun NH2-terminal kinase (1). MEK4 is part of a signaling system that is stimulated in the presence of stress to the cell (including chemotherapy), and it controls differentiation and programmed cell death (apoptosis). MEK4 is central component of signal transduction pathways induced by mitogens and stresses. MEK4 genetic inactivation has been observed in a subset of pancreatic carcinomas, implicating deregulation of the stress-activated protein kinase pathway in pancreatic carcinogenesis (2).

### References

1. Karl Deacon, K. et al: Characterization of the Mitogen-activated Protein Kinase Kinase 4 (MEK4)/c-Jun NH2-terminal kinase 1 and MKK3/p38 Pathways Regulated by MEK Kinases 2 and 3. J. Biol. Chem.1997; 272(22): 14489-14496.
2. Wei, X. et al: MAP2K4/MEK4 Expression in Pancreatic Cancer. Clinical Cancer Res. 2004; 10, 8516-8520.

### Purity



The purity was determined to be **>95%** by densitometry. Approx. MW **63kDa**.

## MEK4 (MAP2K4), Unactive

Recombinant human protein expressed in E. coli cells

Catalog #	M05-15G
Lot #	K305-1
Purity	>95%
Concentration	0.1 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.**