

## MEK3 (MAP2K3), Unactive

Full-length recombinant human protein expressed in Sf9 cells

Catalog # **M04-14G**

Lot # L196-1

### Product Description

Recombinant full-length human MEK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_002756](#).

### Gene Aliases

MAP2K3, MEK3, MAPKK3, PRKMK3, SKK2

### 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.

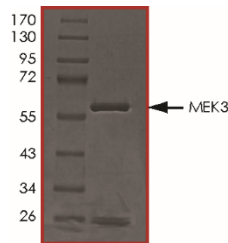
### Scientific Background

Mitogen-activated protein kinase kinase 3 (MEK3) is a serine/threonine kinase that is involved in the activation of MAP kinases. MEK3 is a dual-specificity protein kinase that phosphorylates p38 MAP kinase at its activation site Thr-Gly-Tyr and this stimulates the ability of p38 MAPK to phosphorylate protein substrates such as ATF-2 and Elk-1. MEK3 is activated by different forms of cellular stress and inflammatory cytokines. MEK3-null mice are viable and fertile, but their macrophages and dendritic cells are defective in IL-12 production (1). A virulence factor from *Yersinia pseudotuberculosis*, YopJ, was found to bind directly to MEK3 *in vitro* and this leads to blockage of its phosphorylation and activation (2).

### References

1. Lu, H.-T. et al: Defective IL-12 production in mitogen-activated protein (MAP) kinase kinase 3 (MEK3)-deficient mice. *EMBO J.* 18: 1845-1857, 1999.
2. Orth, K. et al: Inhibition of the mitogen-activated protein kinase kinase superfamily by a *Yersinia* effector. *Science* 285: 1920-1923, 1999.

### Purity



The purity of MEK3 protein was determined to be **>80%** by densitometry, approx. MW **62kDa**.

## MEK3 (MAP2K3), Unactive

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	M04-14G
Lot #	L196-1
Purity	>80%
Concentration	0.1 µg/µl
Stability	1 yr 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.**