

MEK1, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog # M02-14BG

Lot # R038-2

Product Description

Recombinant full-length human MEK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_002755](#).

Gene Aliases

MAP2K1; MKK1; MAPKK1; PRKMK1

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.

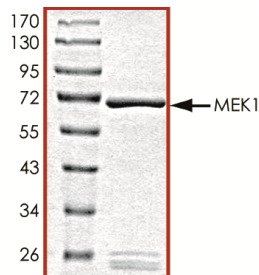
Scientific Background

MEK1 is a protein serine/threonine kinase that is an upstream activator of the extracellular signal-regulated kinases (ERKs) which are activated in response to numerous growth factors and cytokines (1). Activation of ERKs by MEK1 involves phosphorylation of both tyrosine and threonine residues. MEK1, like the ERKs is ubiquitously distributed in tissues with the highest expression in heart, brain and spinal cord. MEK plays an important role in the growth and survival of cancer cells. In hepatocellular carcinoma, MEK activity is frequently elevated and inhibitors of MEK activity may be an effective treatment for this disease (2).

References

1. Sugden, P H. et al: Regulation of the ERK subgroup of MAP kinase cascades through G protein-coupled receptors. *Cell Signal.* 1997 Aug; 9(5):337-51.
2. Huynh, H. et al: Over-expression of the mitogen-activated protein kinase (MAPK) kinase (MEK)-MAPK in hepatocellular carcinoma: its role in tumor progression and apoptosis. *BMC Gastroenterol.* 2003 Aug 8; 3:19.

Purity



The purity of MEK1 was determined to be **>95%** by densitometry. Approx. MW **71kDa**.

Upstream Active Kinase

Active RAF1 (EE)	Cat #R01-13G
Active BRAF	Cat #B08-11G

MEK1, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	M02-14BG
Specific Lot Number	R038-2
Purity	>95%
Concentration	0.2 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
www.signalchem.com

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