

ERK4 (MAPK4) Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # M30-34G

Lot # P1592-4

Product Description

Full-length recombinant human ERK4 (MAPK4) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number for ERK4 (MAPK4) is [NM_002747](#).

Gene Aliases

MAPK4; ERK3; p63MAPK; PRKM4

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.

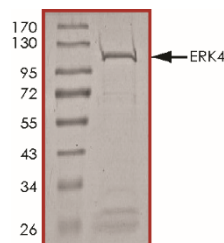
Scientific Background

Extracellular signal-regulated kinase 4 (ERK4) is a serine/threonine protein kinase and belongs to atypical MAPKs. Among MAPKs, it is most closely related to the MAP kinase ERK3, with 73% amino acid identity within the kinase catalytic domain. The precise role of ERK4 is still unclear. It phosphorylates microtubule-associated protein 2 (MAP2) and MAPKAPK5, and may promote entry in the cell cycle.

References

- Rousseau J, et al. Targeted inactivation of Mapk4 in mice reveals specific nonredundant functions of Erk3/Erk4 subfamily mitogen-activated protein kinases. *Mol Cell Biol.* 30:5752-63, 2010
- Kostenko S, et al. Tumour promoting and suppressing roles of the atypical MAP kinase signalling pathway ERK3/4-MK5. *J Mol Signal.* 7:9, 2012.

Purity



The purity of ERK4 (MAPK4) was determined to be **>85%** by densitometry, ERK4 (MAPK4) approx. MW **110kDa**.

ERK4 (MAPK4) Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog#	M30-34G
Lot #	P1592-4
Purity	>85%
Concentration	0.05 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
www.signalchem.com

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