

ERK3 (MAPK6) Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # M31-34G

Lot # A1494-1

Product Description

Recombinant full-length human ERK3 (MAPK6) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM_002748](#).

Gene Aliases

ERK3; MAPK6; HsT17250; p97MAPK; PRKM6

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.

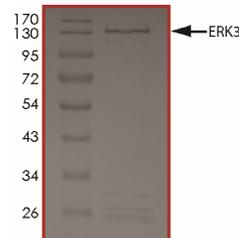
Scientific Background

ERK3 (MAPK6) or mitogen-activated protein kinase 6 is a member of the Ser/Thr protein kinase family, and is most closely related to mitogen-activated protein kinases (MAP kinases) which also known as extracellular signal-regulated kinases (ERKs) that are activated through protein phosphorylation cascades and act as integration points for multiple biochemical signals and tyrosine phosphorylated in response to insulin and NGF (1). ERK3 is highly expressed in various human tissues, most abundantly in skeletal muscle (2).

References

1. Boulton, T. G.et.al: ERKs: a family of protein-serine/threonine kinases that are activated and tyrosine phosphorylated in response to insulin and NGF. Cell 65: 663-675, 1991.
2. Meloche, S. et.al: Primary structure, expression and chromosomal locus of a human homolog of rat ERK3. Oncogene 13: 1575-1579, 1996.

Purity



The purity of ERK3 (MAPK6) was determined to be **>75%** by densitometry. Approx. MW **135 kDa**.

ERK3 (MAPK6) Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	M31-34G
Lot #	A1494-1
Purity	>75%
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.