

MAPKAPK3, Unactive

Full-length recombinant protein expressed in E. coli cells

Catalog # M41-14G

Lot # N200-5

Product Description

Recombinant full-length human MAPKAPK3 was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_004635](#).

Gene Aliases

3PK; MAPKAP3

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.

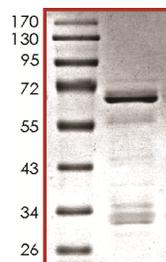
Scientific Background

MAPKAPK3 is closely related to MAPKAPK2 sharing 72% nucleotide and 75% amino acid identity (1). MAPKAPK3 is activated by growth inducers and stress stimulation of cells. *In vitro* studies demonstrated that ERK, p38 MAP kinase and Jun N-terminal kinase were all able to phosphorylate and activate this kinase, which suggested the role of MAPKAPK3 as an integrative element of signaling in both mitogen and stress responses (2).

References

1. Sithanandam, G. et al: 3pK a new mitogen-activated protein kinase-activated protein kinase located in the small cell lung cancer tumor suppressor gene region. *Molec. Cell. Biol.* 16: 868-876, 1996.
2. Ludwig, S. et al: 3pK, a novel mitogen-activated protein (MAP) kinase-activated protein kinase, is targeted by three MAP kinase pathways. *Molec. Cell. Biol.* 16: 6687-6697, 1996.

Purity



← MAPKAPK3 The purity of MAPKAPK3 was determined to be **>90%** by densitometry. Approx. MW **69kDa**.

Upstream Active Kinases

Active P38α	Cat # M39-10
Active P38β	Cat # M36-10
Active P38δ	Cat # M38-10
Active P38γ	Cat # M37-10

MAPKAPK3, Unactive

Full-length recombinant protein expressed in E. coli cells

Catalog Number	M41-14G
Specific Lot Number	N200-5
Purity	>90%
Concentration	0.2 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.