

TOPK, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog # T14-14G

Lot # N319-1

Product Description

Recombinant full-length human TOPK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_018492](#).

Gene Aliases

PBK, SPK, Nori-3, FLJ14385

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.

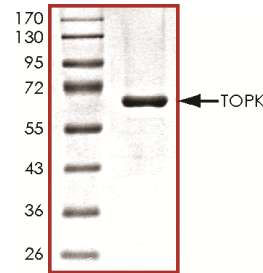
Scientific Background

TOPK is a MAPK kinase that phosphorylates p38 MAPK and is activated in a cell-cycle-dependent manner in neuronal progenitor cells *in vitro* (1). Expression of TOPK is detected in male germ line progenitor cells, activated T-cells, and a variety of lymphomas and leukemias. *In vitro* studies have shown that activated TOPK phosphorylated p38 MAPK but not JNK or ERK. TOPK activation requires phosphorylation by both the M-phase CDK1/CyclinB kinase complex and another unknown kinase, possibly Raf C or Raf A. TOPK may play an important role in linking extracellular signals to an intracellular state, possibly allowing extracellular influence on the cell-cycle-related processes of proliferation or differentiation (2).

References

1. Matsumoto, S. et al: Characterization of a MAPKK-like protein kinase TOPK. *Biochem Biophys Res Commun.* 2004; 325: 997-1004.
2. Simons-Evelyn, M. et al: PBK/TOPK is a novel mitotic kinase which is upregulated in Burkitt's lymphoma and other highly proliferative malignant cells. *Blood Cells Mol Dis.* 2001; 27: 825-829.

Purity



The purity was determined to be **>95%** by densitometry. Approx. MW **68kDa**.

TOPK, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog Number T14-14G

Specific Lot Number N319-1

Purity >95%

Concentration 0.1 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.