

TRIB3 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog # T32-30G

Lot # Y768-3

Product Description

Recombinant full-length human TRIB3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The TRIB3 gene accession number is [NM_021158](#).

Gene Aliases

NIPK, SINK, SKIP3, TRB3

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.

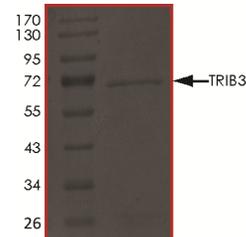
Scientific Background

TRIB3 or tribbles homolog 3 protein encode one of three members of the Tribbles family that share a Trb domain, which is homologous to protein serine-threonine kinases, but lacks the active site lysine and probably lacks a catalytic function. TRIB3 is a negative regulator of NF-kappaB and can also sensitize cells to TNF- and TRAIL-induced apoptosis. TRIB3 can negatively regulate the cell survival serine-threonine kinase AKT1. TRIB3 modulation by H. pylori and/or its products may be an important mechanism of H. pylori-associated pathogenesis (1). TRIB3 is a p65-interacting negative regulator of NF-kappa-B-dependent transcription (2). TRIB3 is highly expressed in pancreas, peripheral blood leukocytes, and bone marrow.

References

1. Qi, L.et.al: TRB3 links the E3 ubiquitin ligase COP1 to lipid metabolism. Science 312: 1763-1766, 2006.
2. Wu, M. et.al: SINK is a p65-interacting negative regulator of NF-kappa-B-dependent transcription. J. Biol. Chem. 278: 27072-27079, 2003.

Purity



The purity of TRIB3 was determined to be **>90%** by densitometry, approx. MW **69kDa**.

TRIB3 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	T32-30G
Lot #	Y768-3
Purity	>90%
Concentration	0.1 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.