

TRIB2 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog # T31-30G

Lot # Y874-1

Product Description

Recombinant full-length human TRIB2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The TRIB1 gene accession number is [NM_021643](#).

Gene Aliases

GS3955, TRB2

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.

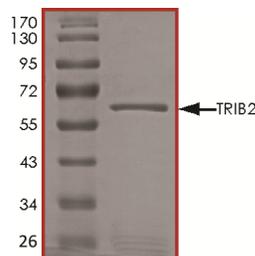
Scientific Background

TRIB2 or tribbles homolog 2 proteins 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. TRIB2 protein interacts and modulates the activity of signal transduction pathways in a number of physiological and pathological processes. TRIB2 is regulated by inflammatory stimuli in myeloid (THP-1) cells. TRIB2 can act as an oncogene that inactivates the transcription factor C/EBPalpha (CCAAT/enhancer-binding protein alpha) and causes acute myelogenous leukemia (1). TRIB2 is a p65-interacting negative regulator of NF-kappa-B-dependent transcription (2).

References

1. Wilkin, F. et.al: Characterization of a phosphoprotein whose mRNA is regulated by the mitogenic pathways in dog thyroid cells. *Europ. J. Biochem.* 248: 660-668, 1997.
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 TRIB2 was determined to be **>90%** by densitometry, approx. MW **64kDa**.

TRIB2 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	T31-30G
Lot #	Y874-1
Purity	>90%
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.