

TBCKL Protein

Full-length recombinant human protein expressed in sf9 cells

Catalog # T35-31G

Lot # U1511-3

Product Description

Full-length recombinant human TBCKL was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC009208](#).

Gene Aliases

HSPC302

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.

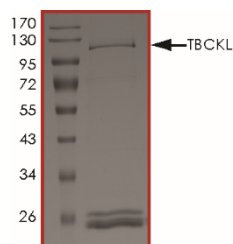
Scientific Background

TBC domain-containing protein kinase-like protein (TBCK) is a Ser/Thr kinase that contains a protein kinase domain, a Rhodanase-like domain and the Tre-2/Bub2/Cdc16 (TBC) domain. TBCK is involved in the regulation of mTOR signaling pathway and plays a role in cell proliferation, cell size, and the organization of actin. 3 isoforms of the human protein are produced by alternative splicing.

References

1. Liu Y, et al: TBCK influences cell proliferation, cell size and mTOR signaling pathway. PLoS One. 2013 Aug 19;8(8):e71349.
2. <http://www.phosphosite.org/proteinAction.do?id=2231&showAllSites=true>

Purity



The purity of TBCKL protein was determined to be **>70%** by densitometry, TBCKL approx. MW **118kDa**.

TBCKL Protein

Full-length recombinant human protein expressed in sf9 cells

Catalog #	T35-31G
Lot #	U1511-3
Purity	>70%
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.