

TPM3 Protein

Full-length human recombinant protein expressed in E.coli cells

Catalog # T36-30H

Lot # Y973-4

Product Description

Full-length recombinant human TPM3 protein was expressed in E.coli cells using an N-terminal His tag. The gene accession number is [NM_152263](#).

Gene Aliases

APM1; CFTD; hscp30; NEM1; OK/SW-cl.5; TM-5; TM3; TM30; TM30nm; TM5; TPMsk3

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 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.

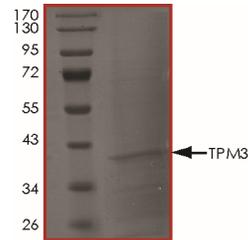
Scientific Background

TPM3 or tropomyosin 3 is a member of the tropomyosin family of actin-binding proteins that are dimers of coiled-coil proteins that provide stability to actin filaments and regulate access of other actin-binding proteins. TPM3 involved in translocations with other loci, including anaplastic lymphoma receptor tyrosine kinase (ALK) and neurotrophic tyrosine kinase receptor type 1 (NTRK1), which result in the formation of fusion proteins that act as oncogenes. TPM3 protein is substituted for the external domain of a putative tyrosine-kinase cell surface receptor to create the TRK oncogene (1). TPM3 are a common cause of congenital fiber type disproportion (2).

References

1. Coulier, F.et.al: Mechanism of activation of the human TRK oncogene. Molec. Cell. Biol. 9: 15-23, 1989.
2. Clarke, N. F. et.al: Mutations in TPM3 are a common cause of congenital fiber type disproportion. Ann. Neurol. 63: 329-337, 2008.

Purity



The purity of TPM3 protein was determined to be **>70%** by densitometry. Approx. MW **39kDa**.

TPM3 Protein

Full-length human recombinant protein expressed in E.coli cells

Catalog #	T36-30H
Lot #	Y973-4
Purity	>70%
Concentration	0.1µg/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.