

TRIM32 Protein

Recombinant protein expressed in Sf9 cells

Catalog # T291-30G

Lot # J715-2

Product Description

Recombinant human TRIM32 (2-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_012210](#).

Gene Aliases

BBS11; HT2A; LGMD2H; TATIP

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

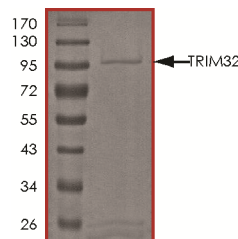
Scientific Background

TRIM32 or tripartite motif containing 32 is a member of the tripartite motif (TRIM) family whose members are involved in diverse cellular functions such as developmental patterning and oncogenesis. TRIM32 localizes to cytoplasmic bodies and nucleus, where it interacts with the activation domain of the HIV-1 Tat protein that activates transcription of HIV-1 genes. TRIM32 is a regulator of dysbindin and mutations in TRIM32 may impair substrate ubiquitination (1). Deficiency of TRIM32 may involve both neurogenic and myogenic characteristics (2).

References

1. Locke, M.et.al: TRIM32 is an E3 ubiquitin ligase for dysbindin. Hum. Molec. Genet. 18: 2344-2358, 2009.
2. Kudryashova, E.et.al: Deficiency of the E3 ubiquitin ligase TRIM32 in mice leads to a myopathy with a neurogenic component. Hum. Molec. Genet. 18: 1353-1367, 2009.

Purity



The purity of TRIM32 was determined to be **>75%** by densitometry. Approx. MW **100 kDa**.

TRIM32 Protein

Recombinant protein expressed in Sf9 cells

Catalog Number	T291-30G
Specific Lot Number	J715-2
Purity	>75%
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.