

TDP43 Protein

Recombinant human protein expressed in E. coli cells

Catalog # T74-30G

Lot # U1536-5

Product Description

Recombinant human TDP43 protein was expressed in E. coli cells using an N-terminal GST tag. The gene accession number for TDP43 is [NM_007375](#).

Gene Aliases

ALS10; TARDBP

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.

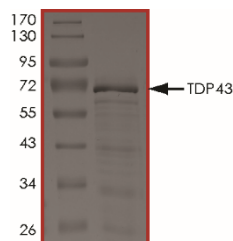
Scientific Background

TAR DNA binding protein (TARDBP) Human was originally discovered in a screen for transcriptional repressors of the transactive response (TAR) DNA binding element of the HIV-1 virus. Contain two RNA-recognition motifs (RRM1 and RRM2), one of which partially overlaps with a putative nuclear export sequence. Mutations in TARDBP were identified in amyotrophic lateral sclerosis (ALS) patients.

References

1. Ou SH, et al: Cloning and characterization of a novel cellular protein, TDP-43, that binds to human immunodeficiency virus type 1 TAR DNA sequence motifs. *J Virol* 69:3584–3596. 1995
2. Gitcho MA, et al. TDP-43 A315T mutation in familial motor neuron disease. *Ann Neurol* 2008 63:535–538. 2008

Purity



The purity of TDP43 protein was determined to be **>70%** by densitometry, TDP43 approx. MW **72kDa**.

TDP43 Protein

Recombinant human protein expressed in E. coli cells

Catalog #	T74-30G
Lot #	U1536-5
Purity	>70%
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.