

Recombinant TRIT1 (48-467) protein

Catalog No: 81190

Expressed In: Baculovirus

Quantity: 20 µg

Concentration: 0.15 µg/µl

Source: Human

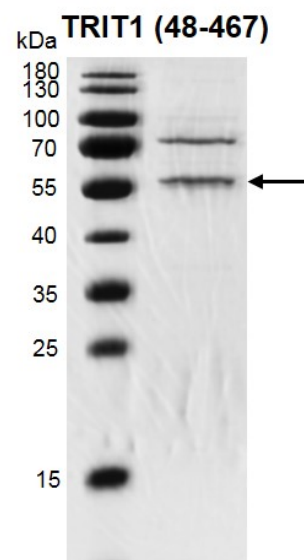
Buffer Contents: Recombinant TRIT1 (48-467) protein is supplied in 25 mM HEPES/NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

Background: TRIT1 (tRNA Isopentenyltransferase 1), also called as IPTase or MOD5, is a highly conserved tRNA isopentenyl transferase that modifies a subset of tRNAs in human cells. It is targeted to the mitochondrion and modifies transfer RNAs (tRNAs) by adding a dimethylallyl group onto the adenine at position 37. This modification is important for maintaining the correct reading frame during protein translation. TRIT1 is considered a tumor suppressor and its expression can decrease cell growth. Alternative splicing results in multiple transcripts variants, most of which are likely non-functional.

Protein Details: Recombinant TRIT1 (48-467) protein that includes amino acids 48-467 of human TRIT1 protein (accession number NP_060116.2) was expressed in a baculovirus expression system and contains an N-terminal FLAG-Tag with a molecular weight of 49.3 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data this product is shown.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



Recombinant TRIT (48-467) protein gel
10% SDS-PAGE gel with Coomassie blue staining
MW: 49.3 kDa
Purity: >55%