

Recombinant WTAP protein

Catalog No: 31571, 31971

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.2 µg/µl

Source: Human

Buffer Contents: Recombinant WTAP protein is supplied at a concentration of 0.2 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

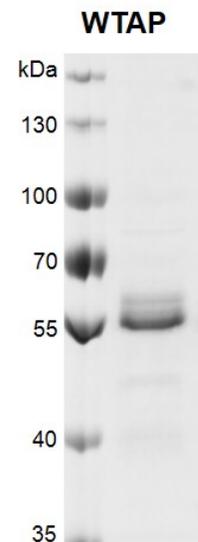
Background: N6-methylated adenine (m6A) is prevalently present in nearly all RNA types and can be found in all organisms from bacteria to humans. It preferentially appears around stop codons and within long internal exons in mammalian messenger RNAs. m6A plays an important role in the efficiency of mRNA splicing, processing, translation efficiency, editing and mRNA stability. m6A also takes place in other RNA molecules, such as primary miRNA (pri-miRNAs).

WTAP (Wilms Tumor 1 Associated Protein, as known as Mum2) is a ubiquitously expressed nuclear protein. WTAP is the regulatory subunit of the WMM N6-methyltransferase complex, a multiprotein complex that mediates N6-methyladenosine (m6A) methylation of some adenosine residues of some mRNAs. WTAP is required for accumulation of METTL3 and METTL14 to nuclear speckle. It acts as a mRNA splicing regulator. WTAP also regulates G2/M cell-cycle transition by binding to the 3' UTR of CCNA2, which enhances its stability. WTAP can impair WT1 DNA-binding ability and inhibit expression of WT1 target genes.

Protein Details: Recombinant WTAP protein was expressed in a baculovirus expression system as the full length protein (accession number NP_004897.2) with an N-terminal FLAG tag. The molecular weights of WTAP is 45.2 kDa.

Application Notes: Recombinant WTAP protein is suitable for use in protein-protein interaction, RNA binding assay.

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 WTAP protein gel. Recombinant WTAP protein was run on an 8% SDS-PAGE gel and stained with Coomassie blue.