

Recombinant METTL14 protein

Catalog No: 31568, 31968

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.2 µg/µl

Source: Human

Buffer Contents: Recombinant METTL14 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.

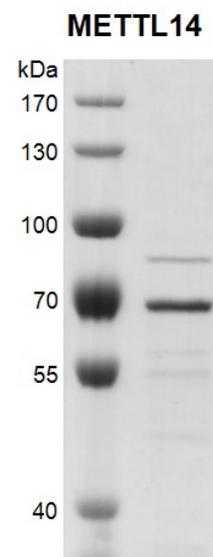
Background: N6-methylated adenine (m6A) is prevalently present in nearly all RNA types and can be found in all organisms from bacteria to humans. It preferentially appears around stop codons and within long internal exons in mammalian messenger RNAs. m6A plays an important role in the efficiency of mRNA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

METTL14 (methyltransferase-like 14) forms a stable N6-methyltransferase heterodimer complex with METTL3, which catalyzes the generation of m6A modification on mammalian nuclear RNAs. METTL3 is the catalytically active subunit, while METTL14 plays a structural role critical for substrate recognition.

Protein Details: Recombinant METTL14 protein was expressed in a Baculovirus expression system as the full length protein (accession number NP_066012.1) with an N-terminal FLAG tag. The molecular weight of METTL14 is 53.3 kDa.

Application Notes: Recombinant METTL14 protein is suitable for use in enzyme kinetics, inhibitor screening, and selectivity profiling.

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 METTL14 protein gel
Recombinant METTL14 was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.

MW: 53.3 kDa

Purity: > 75%