

## Recombinant METTL3 protein

**Catalog No:** 31567, 31967

**Lot No:** 30716001

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.35 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant METTL3 protein is supplied at a concentration of 0.35 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

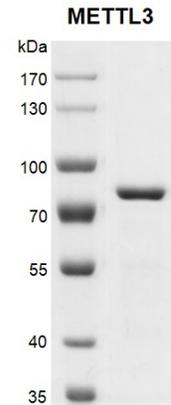
**Background:** N6-methylated adenine (m6A) is prevalently present in nearly all RNA types and can be found in all organisms from bacteria to humans. It preferentially appears around stop codons and within long internal exons in mammalian messenger RNAs. m6A plays an important role in the efficiency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

METTL3 (methyltransferase-like 3, also known as IME4, M6A, MT-A70) forms a stable N6-methyltransferase heterodimer complex with METTL14, 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. METTL3 promotes translation by interacting with the translation initiation machinery in the cytoplasm. Its overexpression in a number of cancer cells suggests that it may participate to cancer cell proliferation by promoting mRNA translation.

**Protein Details:** Recombinant METTL3 protein was expressed in a Baculovirus expression system as the full length protein (accession number NP\_062826.2) with an N-terminal FLAG-Tag. The molecular weight of METTL3 is 65.5 kDa.

**Application Notes:** Recombinant METTL3 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 METTL3 protein gel.**

Recombinant METTL3 protein was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.

MW: 65.5 kDa

Purity: > 95%