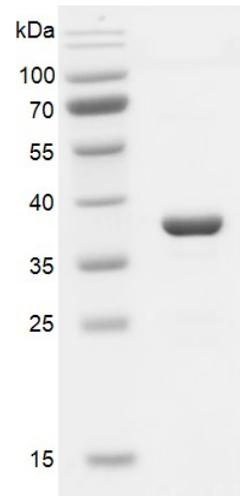


Recombinant METTL1 protein, His-Tag

Catalog No: 81058, 81758**Lot No:** 27017001**Expressed In:** *E. coli***Quantity:** 100, 1000 µg**Concentration:** 1.5 µg/µl**Source:** Human**Buffer Contents:** Recombinant METTL1 protein, His-Tag is supplied at a concentration of 1.5 µg/µl in 25 mM Tris pH 8.0, 300 mM NaCl, 10% glycerol, 0.5 mM TCEP.**Background:** METTL1 (methyltransferase-like 1, also known as TRM8, TRMT8, C12orf1, YDL201w) contains a conserved S-adenosylmethionine-binding motif and is inactivated by phosphorylation. METTL1 forms a stable heterodimer complex with WDR4, which catalyzes the formation of N(7)-methylguanine at position 46 (m7G46) in tRNA.**Protein Details:** Recombinant METTL1 protein, His-Tag was expressed in *E. coli* cells as the full length protein (accession number NP_005362.3) with an N-terminal His-Tag. The molecular weight of METTL1, His-Tag is 35 kDa**Application Notes:** Recombinant METTL1 protein, His-Tag 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.

His-METTL1

**Recombinant METTL1, His-Tag, protein gel.**

12.5% SDS-PAGE gel, stained with Coomassie blue.

MW: 35 kDa

Purity: ≥ 95%