

## Recombinant METTL1 protein

**Catalog No:** 81022

**Lot No:** 14417001

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.2 µg/µl

**Source:** Human

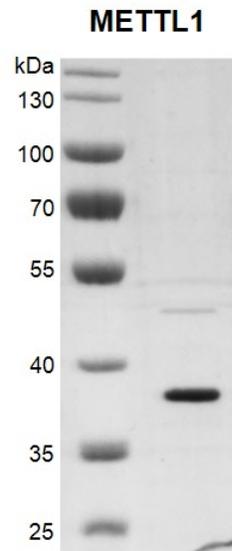
**Buffer Contents:** Recombinant METTL1 protein is supplied at a concentration of 0.2 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

**Background:** 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 was expressed in a baculovirus expression system as the full length protein (accession number NP\_005362.3) with an N-terminal FLAG tag. The molecular weights of METTL1 is 32.7 kDa.

**Application Notes:** Recombinant METTL1 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 METTL1 protein gel**

10% SDS-PAGE Coomassie staining

MW: 32.7 ka

Purity: ≥ 87%