

## Recombinant METTL2A protein

**Catalog No:** 81027, 81727

**Lot No:** 22517001

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.3 µg/µl

**Source:** Human

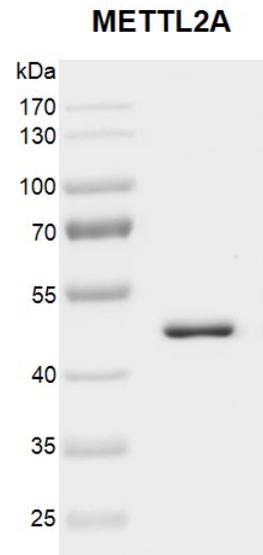
**Buffer Contents:** Recombinant METTL2A protein is supplied at a concentration of 0.3 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

**Background:** METTL2A (methyltransferase-like 2A) contains a conserved S-adenosylmethionine-binding motif. METTL2A is a probable S-adenosyl-L-methionine-dependent methyltransferase that mediates 3-methylcytidine modification of some tRNAs

**Protein Details:** Recombinant METTL2A protein was expressed in a baculovirus expression system as the full length protein (accession number NP\_859076.3) with an N-terminal FLAG tag. The molecular weights of METTL2A is 44.8 kDa.

**Application Notes:** Recombinant METTL2A 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 METTL2A protein gel**  
10% SDS-PAGE Coomassie staining  
MW: 44.8 kDa  
Purity: ≥95%