

Recombinant METTL18 protein

Catalog No: 81030, 81730

Lot No: 23817001

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.3 µg/µl

Source: Human

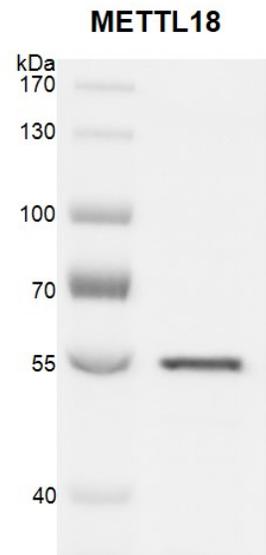
Buffer Contents: Recombinant METTL18 protein is supplied at a concentration of 0.3 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

Background: METTL18 (methyltransferase-like 18, also known as HPM1, AsTP2, C1orf156) contains a conserved S-adenosylmethionine-binding motif. It is a potential SAM-dependent methyltransferase.

Protein Details: Recombinant METTL18 protein was expressed in a baculovirus expression system as the full length protein (accession number NP_219486.1) with an N-terminal FLAG tag. The molecular weights of METTL18 is 43.4 kDa.

Application Notes: Recombinant METTL18 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 METTL18 protein gel
12.5% SDS-PAGE Coomassie staining
MW: 43.4 kDa
Purity: ≥95%