

Recombinant METTL4 protein

Catalog No: 81177, 81877

Expressed In: Baculovirus

Quantity: 10 µg

Concentration: 0.15 µg/µl

Source: Human

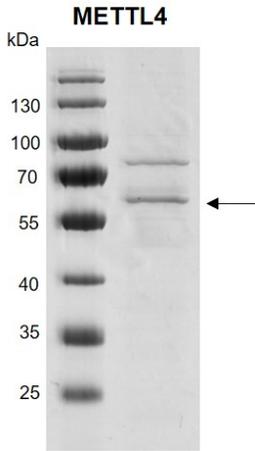
Buffer Contents: Recombinant METTL4 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

Background: METTL4 (Methyltransferase Like 4), is a member of METTL family proteins. Gene Ontology (GO) annotations related to this gene include nucleic acid binding and methyltransferase activity. It is a probable methyltransferase.

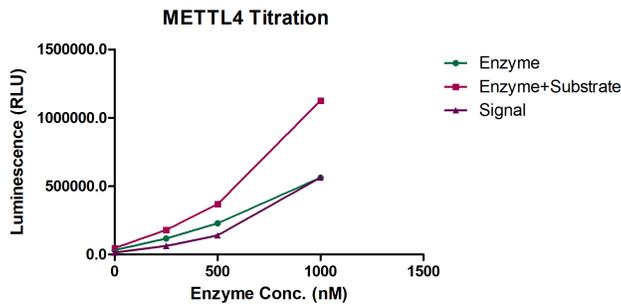
Protein Details: Recombinant METTL4 protein was expressed in baculovirus expression system as the full length protein (accession number NP_073751.3) with an N-terminal FLAG-Tag. The molecular weight of the protein is 55.2 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

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.

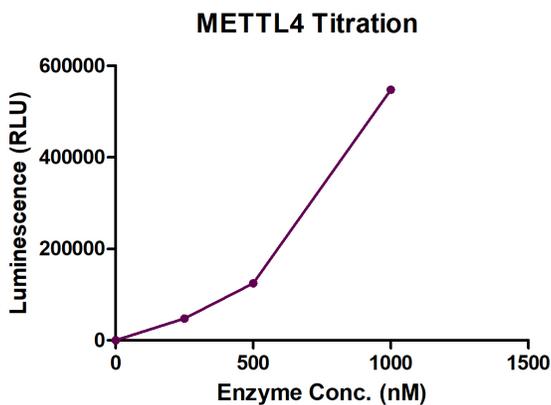


Recombinant METTL4 protein
 10% SDS-PAGE Coomassie staining
 MW: 55.2 kDa
 Purity: >50%



MTase-Glo assay for METTL4 methyltransferase activity

400 ng total RNA and 1 μ M SAM was incubated with different concentrations of METTL4 protein in 8 μ l reaction system containing 50 mM Tris-HCl pH 8.6, 0.02% Triton X -100, 2 mM MgCl₂, 1 mM TCEP at room temperature for 1 hour (0.2 U/ μ L RRI was added in this system). 5 \times MTase-Glo Reagent was added to the products and incubated for 30 min. Then MTase-Glo Detection was added, and luminescence was read after another 30 min incubation.



MTase-Glo assay for METTL4 methyltransferase activity

400 ng total RNA and 1 μ M SAM was incubated with different concentrations of METTL4 protein in 8 μ l reaction system containing 50 mM Tris-HCl pH 8.6, 0.02% Triton X -100, 2 mM MgCl₂, 1 mM TCEP at room temperature for 1 hour (0.2 U/ μ L RRI was added in this system). 5 \times MTase-Glo Reagent was added to the products and incubated for 30 min. Then MTase-Glo Detection was added, and luminescence was read after another 30 min incubation.