

Recombinant METTL8 protein, GST-Tag

Catalog No: 81106, 81806

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 0.5 µg/µl

Source: Human

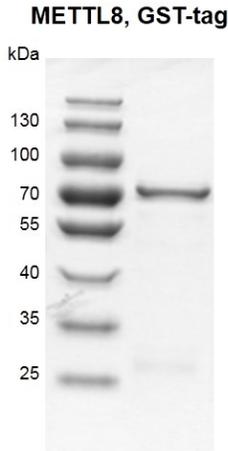
Buffer Contents: Recombinant METTL8 protein, GST-tag protein is supplied in 25 mM Tris-HCl pH 8.0, 500 mM NaCl, 10% glycerol and 0.5 mM TCEP.

Background: METTL8 (methyltransferase-like 8, also known as TIP) contains a conserved S-adenosylmethionine-binding motif. It is a potential SAM-dependent methyltransferase.

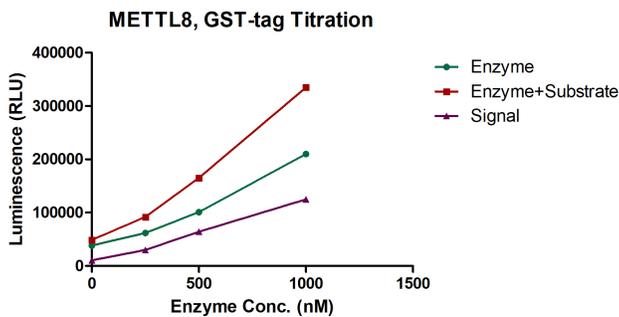
Protein Details: Full length recombinant METTL8 (accession number NP_079046.2) was expressed in *E. coli* cells with an N-terminal GST tag. The molecular weight of the protein is 73.3 kDa.

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.

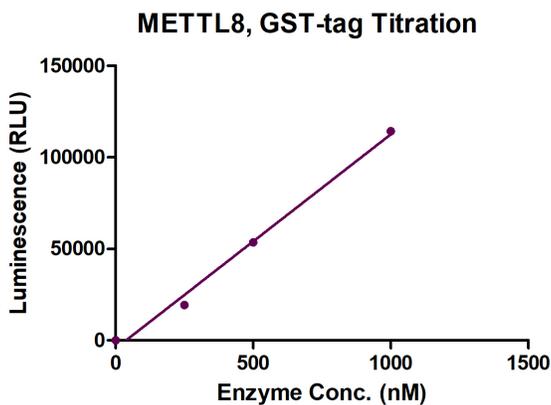


Recombinant METTL8 protein, GST-tag
 10% SDS-PAGE Coomassie staining
 MW: 73.3 kDa
 Purity: ≥90%



MTase-Glo assay for METTL8 methyltransferase activity

400 ng total RNA and 1 μ M SAM was incubated with different concentrations of METTL8 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 METTL8 methyltransferase activity

400 ng total RNA and 1 μ M SAM was incubated with different concentrations of METTL8 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.