

## Recombinant METTL13 protein

---

**Catalog No:** 81024, 81724

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.4 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant METTL13 protein is supplied at a concentration of 0.4 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

**Background:** METTL13 (methyltransferase-like 13, is known as feat, CGI-01, KIAA0859, 5630401D24Rik) contains a conserved S-adenosylmethionine-binding motif, It is a potential SAM-dependent methyltransferase.

**Protein Details:** Recombinant METTL13 protein was expressed in a baculovirus expression system as the full length protein (accession number NP\_057019.3) with an N-terminal FLAG tag. The molecular weights of METTL13 is 80 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 -20°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.

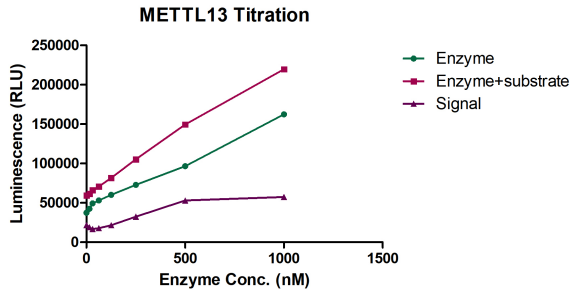


**Recombinant METTL13 protein gel.**

10% SDS-PAGE Coomassie staining

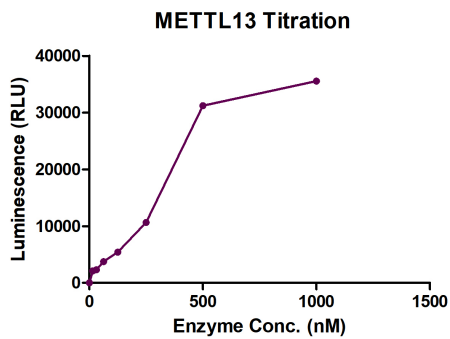
MW: 80 kDa

Purity:  $\geq 87\%$



**MTase-Glo assay for METTL13 methyltransferase activity**

1  $\mu\text{M}$  EEF1A1 protein and 1  $\mu\text{M}$  SAM was incubated with different concentrations of METTL13 protein in 8  $\mu\text{l}$  reaction system containing 50 mM Tris-HCl pH 8.6, 0.02% Triton X-100, 2 mM MgCl<sub>2</sub>, 1 mM TCEP at room temperature for 1 hour. 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 METTL13 methyltransferase activity**

1  $\mu\text{M}$  EEF1A1 protein and 1  $\mu\text{M}$  SAM was incubated with different concentrations of METTL13 protein in 8  $\mu\text{l}$  reaction system containing 50 mM Tris-HCl pH 8.6, 0.02% Triton X-100, 2 mM MgCl<sub>2</sub>, 1 mM TCEP at room temperature for 1 hour. 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.