

Recombinant METTL7B protein

Catalog No: 81208, 81908

Expressed In: Baculovirus

Quantity: 20 µg

Concentration: 0.2 µg/µl

Source: Human

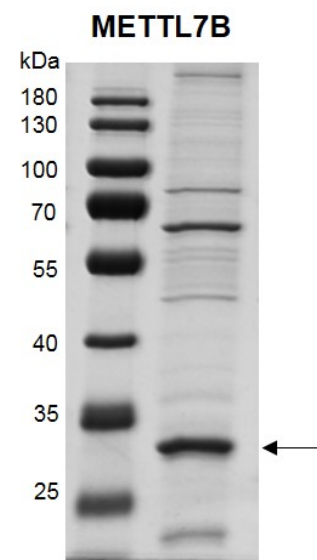
Buffer Contents: Recombinant METTL7B protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

Background: METTL7B (Methyltransferase-Like Protein 7B) was originally identified as a Golgi-associated methyltransferase and is S-adenosylmethionine-dependent. It is associated with a lot of diseases including Muscular Dystrophy, Congenital, Due To Integrin Alpha-7 Deficiency and Supraglottis Cancer. Previous work has shown that RhoBTB1 regulates the integrity of the Golgi complex through METTL7B. RhoBTB1 is required for expression of METTL7B and silencing of either protein leads to fragmentation of the Golgi.

Protein Details: Recombinant METTL7B protein was expressed in a baculovirus expression system as the full length human protein (accession number NP_004681.1) with a C-terminal FLAG tag. The molecular weight of the protein is 29.5 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 this product is shown.

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 METTL7B protein gel
10% SDS-PAGE gel with Coomassie blue staining
MW: 29.5 kDa
Purity: >65%