

Recombinant METTL22 protein

Catalog No: 81092, 81892

Lot No: 31617001

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.2 µg/µl

Source: Human

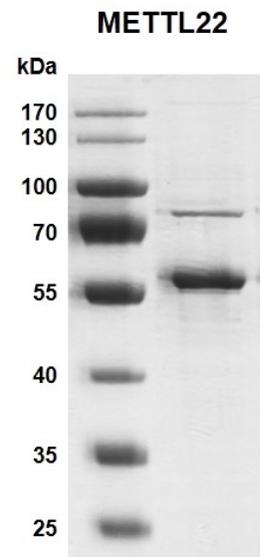
Buffer Contents: Recombinant METTL22 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100 and 0.5 mM TCEP.

Background: METTL22 (Methyltransferase Like 22), also called C16orf68 (Chromosome 16 Open Reading Frame 68) or LP8272, is a member of the non-histone lysine methyltransferases. It maybe a S-adenosylmethionine-dependent methyltransferases. *In vitro*, METTL22 interacts with its substrate, Kin17, which is involved in DNA repair and replication and mRNA processing.

Protein Details: Recombinant human METTL22 protein was expressed in a baculovirus expression system as the full length protein (accession number NP_077014.3) with an N-terminal FLAG tag. The molecular weight of the protein is 45.8 kDa.

Application Notes: This protein is suitable for use in binding assays, 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 METTL22 protein gel
10% SDS-PAGE Coomassie staining
MW: 45.8 kDa
Purity: >80%