

Recombinant METTL19 (TRMT44) protein

Catalog No: 81088, 81788

Lot No: 31617001

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.35 µg/µl

Source: Human

Buffer Contents: Recombinant METTL19 (TRMT44) 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.

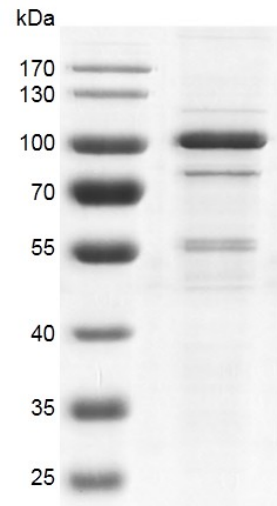
Background: METTL19 (Methyltransferase-Like Protein 19), also known as TRMT44 (TRNA Methyltransferase 44 Homolog), is a putative tRNA methyltransferase found in the cytoplasm. It has a probable adenosyl-L-methionine (AdoMet)-dependent tRNA (uracil-O(2)-)-methyltransferase activity. Defects in *METTL19* gene may be associated with partial epilepsy with pericentral spikes (PEPS).

Protein Details: Recombinant human METTL19 (TRMT44) protein was expressed in a baculovirus expression system as the full length protein (accession number NP_689757.2) with an N-terminal FLAG tag. The molecular weight of the protein is 85.8 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 guaranteed for 6 months from date of arrival.

METTL19 (TRMT44)



Recombinant METTL19 (TRMT44) protein gel
10% SDS-PAGE Coomassie staining
MW: 85.8 kDa
Purity: >75%