

METTL14 Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog # M334-30G

Lot # B2033-7

Product Description

Recombinant full-length human METTL14 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_020961](#).

Gene Aliases

(None)

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

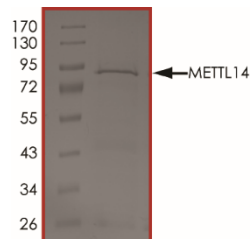
Scientific Background

N6-adenosine-methyltransferase subunit METTL14 (Methyltransferase-like protein 14 or METTL14) methylates adenosine residues of some mRNAs and acts as a regulator of the circadian clock and differentiation of embryonic stem cells. N6-methyladenosine (m6A), which takes place at the 5'-[AG] GAC-3' consensus sites of some mRNAs, plays a role in the efficiency of mRNA splicing, processing and mRNA stability.

References

1. <http://www.uniprot.org/uniprot/Q9HCE5>
2. Liu J, et al: METTL3-METTL14 complex mediates mammalian nuclear RNA N6-adenosine methylation. Nat Chem Biol. 10(2):93-5, 2014.

Purity



The purity of METTL14 was determined to be **>85%** by densitometry. Approx. MW **~84kDa**.

METTL14 Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog Number	M334-30G
Lot #	B2033-7
Purity	>85%
Concentration	0.05 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.