

## METTL3 Protein

Full length human recombinant protein expressed in Sf9 cells

### Catalog # M323-30G

Lot # B1997-6

### Product Description

Recombinant full-length human METTL3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_019852](#).

### Gene Aliases

IME4; M6A; MT-A70; Spo8

### 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.

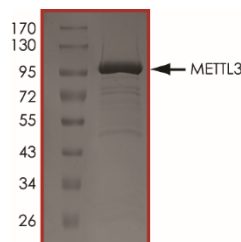
### Scientific Background

N6-adenosine-methyltransferase 70 kDa subunit (MT-A70 or METTL3) belongs to the methyltransferase superfamily that methylates adenosine residues of some RNAs. It acts as a regulator of the circadian clock, differentiation of embryonic stem cells and primary miRNA processing.

### References

1. <http://www.uniprot.org/uniprot/Q86U44>
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 METTL3 was determined to be **>95%** by densitometry. Approx. MW **~100kDa**.

## METTL3 Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog Number	M323-30G
Lot #	B1997-6
Purity	>95%
Concentration	0.1 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.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**