

METTL3/METTL14, Active

Recombinant human proteins expressed in Sf9 cells

Catalog # M323-380G

Lot # B2132-12

Product Description

Full-length recombinant human METTL3/METTL14 complex was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The METTL3 and METTL14 gene accession numbers are [NM_019852](#) and [NM_020961](#).

Gene Aliases

METTL3: IME4; M6A; MT-A70; Spo8

METTL14: None

Formulation

Recombinant proteins 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.

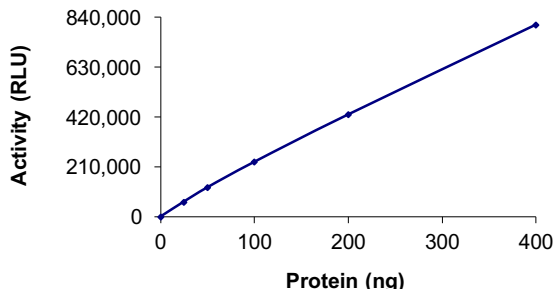
Scientific Background

Methyltransferase-like 3 (METTL3) and methyltransferase-like 14 (METTL14) form a stable heterodimer core complex that catalyzes N⁶-methyladenosine (m⁶A) RNA methylation in mammalian cells. m⁶A is an abundant internal modification in messenger RNA and long non-coding RNA. It functions in multiple aspects of developmental regulation, cell cycle progression, cell fate, and the heat shock stress response by affecting aspects of RNA metabolism such as pre-mRNA processing, translation efficiency, transcript stability and miRNA biogenesis.

References

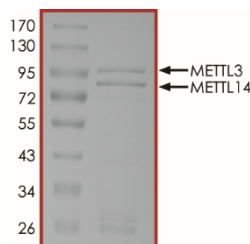
1. Liu J. et al: A METTL3–METTL14 complex mediates mammalian nuclear RNA N⁶-adenosine methylation. *Nat. Chem. Biol.* 2014; 10: 93–95
2. Wang X. et al: Structural basis of N⁶-adenosine methylation by the METTL3–METTL14 complex. *Nature* 2016; 534:575-578
3. Wang P. et al: Structural Basis for Cooperative Function of Mettl3 and Mettl14 Methyltransferases. *Mol. Cell* 2016; 63(2):306-317

Specific Activity



The specific activity of METTL3/METTL14 was determined to be **750 pmol /min/mg** as per activity assay protocol.

Purity



The purity of METTL3/METTL14 complex was determined to be **>90%** by densitometry, approx. MW **100 kDa (METTL3)** and **84 kDa (METTL14)**.

METTL3/METTL14, Active

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	M323-380G
Specific Activity	750 pmol/min/mg
Lot #	B2132-12
Purity	>90%
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
www.signalchem.com

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

Activity Assay Protocol

Reaction Components

Active Methyltransferase (Catalog #: M323-380G)

Active METTL3/METTL14 complex (0.1µg/µl) diluted with RNA MTase Reaction Buffer and assayed as outlined in sample activity plot. (Note: these are suggested working dilutions and it is recommended that the researcher perform a serial dilution of Active MTase complex for optimal results).

RNA Methyltransferase Reaction Buffer

Buffer components: 20mM Tris-HCl, pH7.5, 0.01% Triton X100. Add 1mM DTT (Signalchem, Catalog #: D86-09B-10) and 1unit/µl RNasin® Plus (Promega, Catalog #: N2611, optional) prior to use.

MTase-Glo™ Methyltransferase Assay (Promega, Catalog #: V7601)

S-Adenosyl-Methionine (SAM), 1mM
S-Adenosyl-Homocysteine (SAH), 15 µM
Methyltransferase-Glo™ Reagent, 10X
MTase-Glo™ Detection Solution, 1 bottle

Substrate (Catalog #: M323-58)

METTL3/METTL14 Substrate, an oligo ssRNA, was reconstituted in 25µl of Methyltransferase Dilution Buffer IV to a final concentration of 100µM.

Assay Protocol

The METTL3/METTL14 assay is performed using the Methyltransferase-Glo™ Assays kit (Promega).

- Step 1.** Thaw the active MTase complex and all Methyltransferase-Glo™ Assays kit reagents on ice.
- Step 2.** Prepare the following working solutions with RNA Methyltransferase Reaction Buffer:
 - o 2X final concentration of Active complex (Catalog # M323-380G)
 - o 2X Substrate Cocktail: 40µM of SAM + 4µM of METTL3/METTL14 Substrate (Catalog #: M323-58)
- Step 3.** In a polystyrene 96-well plate, add the following components to bring the initial reaction volume to 20 µl:
 - Component 1.** 10 µl of 2X Substrate Cocktail
 - Component 2.** 10 µl of 2X Active METTL3/METTL14 complex

Note: A blank control can be set up as outlined in step 3 by replacing the enzyme working solution with an equal volume of RNA Methyltransferase Reaction Buffer.

- Step 4.** Mix the reaction on an orbital shaker for 2 minutes. Seal the plate with a plate seal and incubate at 37°C for 60 minutes
- Step 5.** Dilute 10X Methyltransferase-Glo™ Reagent with equal volume of nanopure water, and add 5 µl of the 5X Methyltransferase-Glo™ Reagent to all reaction wells
- Step 6.** Mix on an orbital shaker for 2 minutes and then incubate at room temperature for 30 minutes.
- Step 7.** Add 25 µl of MTase-Glo™ Detection Solution to all reaction wells. Mix for 2 minutes and then incubate at room temperature for 30 minutes
- Step 8.** Read the plate using the KinaseGlo Luminescence Protocol on a GloMax plate reader (Promega; Cat# E7031)
- Step 9.** Using the SAH standard curve, determine the concentration of SAH produced (nM) and calculate the methyltransferase specific activity as outlined below. For a detailed protocol of how to determine SAH amount from RLUs, see MTase-Glo™ Methyltransferase Assay protocol at Promega's website: www.promega.com/protocols

Methyltransferase Specific Activity (SA) (nmol/min/mg)

$$= \frac{[SAH](nM) \times Reaction Volume(\mu l)}{Reaction Time (min) \times Enzyme Amount (mg)} \times 10^{-6}$$

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

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