

## LCMT2 Protein

Recombinant full-length human protein expressed in Sf9 cells

**Catalog # L322-30G**

Lot # B2241-2

### Product Description

Recombinant full-length human LCMT2 protein was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The LCMT2 gene accession number is [NM\\_014793](#).

### Gene Aliases

PPM2; TYW4

### 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.

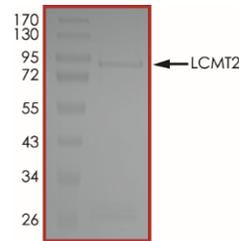
### Scientific Background

LCMT2, also named tRNA wybutosine-synthesizing protein 4 (TYW4), is an S-adenosylmethionine (SAM)-dependent enzyme that involves in the final step of wybutosine biosynthesis pathway (1). Structural study revealed that TYW4 uses one catalytic site for both methylation and methoxycarbonylation process (2).

### References

1. Noma, A., et al. Biosynthesis of wybutosine, a hyper-modified nucleoside in eukaryotic phenylalanine tRNA. EMBO J. 25(10):2142-54, 2006.
2. Suzuki, Y., et al. Structural basis of tRNA modification with CO<sub>2</sub> fixation and methylation by wybutosine synthesizing enzyme TYW4. Nucleic Acids Res. 37(9):2910-25, 2009.

### Purity



The purity of LCMT2 was determined to be **>70%** by densitometry, approx. MW **90 kDa**.

## LCMT2 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	L322-30G
Lot #	B2241-2
Purity	>70%
Concentration	0.05 µg/µl
Stability	1 yr 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.**