

## PRMT7 Protein

Recombinant full-length human protein expressed in Sf9 cells

**Catalog # P365-30GG**

Lot # P1809-3

### Product Description

Recombinant full-length human PRMT7 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The PRMT7 gene accession number is [BC000146](#).

### Gene Aliases

FLJ10640, KIAA1933

### 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.

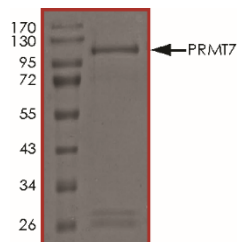
### Scientific Background

PRMT7 (protein arginine methyltransferase 7), an arginine methyltransferase that exhibited type III enzymatic activity, catalyzing the formation of  $\alpha$ -NG-monomethylarginine residues. Specifically mediates the symmetrical dimethylation of arginine residues in the small nuclear ribonucleoproteins Sm D1 (SNRPD1) and Sm D3 (SNRPD3); such methylation being required for the assembly and biogenesis of snRNP core particles. Specifically mediates the symmetric dimethylation of histone H4 'Arg-3' to form H4R3me2s. Plays a role in gene imprinting by being recruited by CTCFL at the H19 imprinted control region (ICR) and methylating histone H4 to form H4R3me2s, possibly leading to recruit DNA methyltransferases at these sites. May also play a role in embryonic stem cell (ESC) pluripotency.

### References

1. Miranda TB, et al: PRMT7 is a member of the protein arginine methyltransferase family with a distinct substrate specificity. *J Biol Chem.* 279(22):22902-7, 2004.
2. Lee JH, et al: PRMT7, a new protein arginine methyltransferase that synthesizes symmetric dimethylarginine. *J Biol Chem.* 280(5):3656-64, 2005.
3. <http://www.uniprot.org/uniprot/Q9NVM4>

### Purity



The purity of PRMT7 was determined to be **>85%** by densitometry, approx. MW **110 kDa**.

## PRMT7 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	P365-30GG
Lot #	P1809-3
Purity	>85%
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.**