

PRMT5 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog # P365-30EG

Lot # Y851-1

Product Description

Recombinant full-length human PRMT5 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The PRMT5 gene accession number is [NM_006109](#).

Gene Aliases

HRMT1L5; IBP72; JBP1; SKB1; SKB1Hs

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.

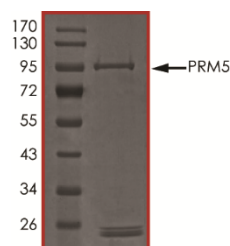
Scientific Background

PRMT5 is a protein arginine methyltransferase and part of a 20S methyltransferase complex that modifies specific arginines to dimethylarginines in several spliceosomal Sm proteins (1). This modification targets Sm proteins to the survival of motor neurons (SMN) complex for assembly into small nuclear ribonucleoprotein core particles. MEP50 associates with PRMT5 in the 20S arginine methyltransferase complex in HeLa cells (2). Antibodies directed against MEP50 can reduce the arginine methyltransferase activity of the immunopurified complex toward Sm substrates. PRMT5-mediated transcriptional repression of reporter gene is dependent on MEP50.

References

1. Friesen, W. J. et al: A novel WD repeat protein component of the methylosome binds Sm proteins. J. Biol. Chem. 277: 8243-8247, 2002.
2. Furuno, K. et al: Association of Polycomb group SUZ12 with WD-repeat protein MEP50 that binds to histone H2A selectively in vitro. Biochem. Biophys. Res. Commun. 345: 1051-1058, 2006.

Purity



The purity of PRMT5 was determined to be **>80%** by densitometry, approx. MW **96 kDa**.

PRMT5 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	P365-30EG
Lot #	Y851-1
Purity	>80%
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.