

PRMT5 recombinant protein

Other names: SKB1, HRMT1L5, SKB1Hs, IBP72, JBP1

Cat. No. C23010107

Source: SF9 cells

Lot #: 001

Size: 10 µg/ 20 µl

Concentration: 0.5 µg/µl

Specificity: Human

Purity: Purified using FPLC, >95% purity as determined by SDS-PAGE.

Storage buffer: 20 mM Tris-Cl pH 7.6, 1 mM EDTA, 0.15 M NaCl, 10% glycerol, 0.5 mM PMSF and 1 mM DTT.

Storage: Store at -80°C; guaranteed stable for 2 years from date of receipt when stored properly.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Description: Full length recombinant human N-terminal Flag tagged PRMT5 (Protein Arginine Methyltransferase 5) protein, produced in SF9 insect cells.

Protein description

PRMT5 (UniProt/Swiss-Prot entry O14744) is an arginine methyltransferase that can both catalyze the monomethylation and symmetrical dimethylation of arginine residues with a preference for the formation of monomethylarginine, in a variety of proteins including SNRPD1, SNRPD3, SUPT5H, MBP, PIWIL1, PIWIL2, PIWIL4, HOXA9, SRGAP2 and histone H2A, H4 and H3. Histone H2A and H4 are methylated at R3 during germ cell development, whereas histone H3 is methylated at R8, probably leading to transcriptional repression.

Quality control

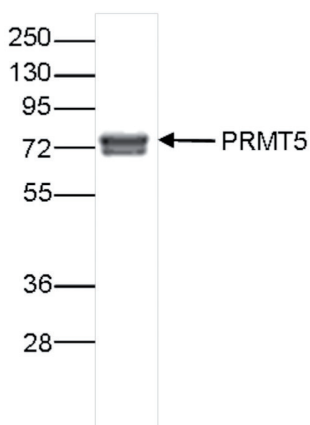


Figure 1.

SDS page of the PRMT5 recombinant protein. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: August 25, 2015