

## PRMT6 (His)

## (Protein Arginine Methyltransferase 6)

**CATALOG NO.:** HMT-21-380

**LOT NO.:**

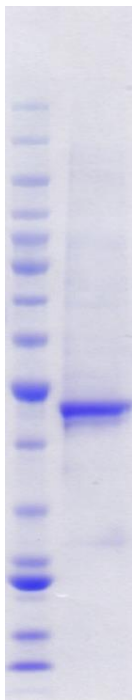
**DESCRIPTION:** Full-length human recombinant PRMT6 expressed in baculovirus system (residues 2-375); Genbank Accession # NM\_018137; N-terminal His-tag; MW = 45.1 kDa). PRMT6, a type I arginine methyltransferase, catalyzes the transfer of a methyl group from S-adenosyl-L-methionine (SAM) to an  $\omega$ -nitrogen of the guanidino function of protein L-arginine residues and the transfer of a second methyl group to the same nitrogen, yielding asymmetric dimethylarginine (aDMA). PRMT6 methylates Arg-2 of histone H3 (H3R2), promoting transcriptional repression.

**PURITY:** >85% by SDS-PAGE.

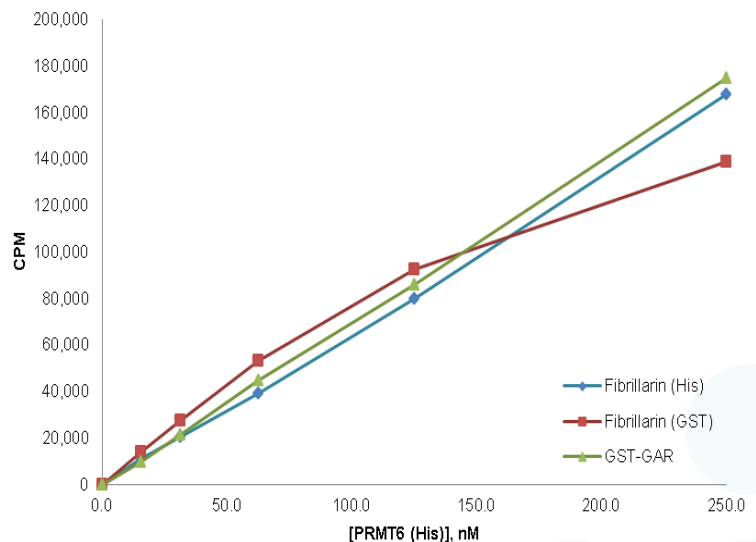
**ASSAY CONDITIONS:** RBC's PRMT6 displays substantial methyltransferase activity from [<sup>3</sup>H]-SAM to GST-GAR (Cat. #HMT-11-137), Fibrillarlin (Hist) (Cat. #HMT-11-183) and Fibrillarlin (GST) (Cat #HMT-11-184; see Figure, below). Activity was determined as TCA-precipitated counts in a scintillation/filter plate assay (Multiscreen FB, Topcount). Reaction conditions: 50 mM Tris-HCl, pH 8.5, 50 mM NaCl, 5 mM MgCl<sub>2</sub>, 1 mM DTT, 1 mM PMSF, 30°C, 60 min. with 1  $\mu$ M [<sup>3</sup>H]-SAM and 1  $\mu$ M of indicated substrate.

**SUPPLIED AS:** \_\_\_  $\mu$ g/ $\mu$ l total protein in 25.4 mM Na<sub>2</sub>HPO<sub>4</sub>, 4.4 mM KH<sub>2</sub>PO<sub>4</sub>, pH 7.4, 137 mM NaCl, 2.7 mM KCl, 3 mM DTT, 30% (w/v) glycerol as determined by OD<sub>280</sub>

**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted enzyme should be refrozen quickly by, for example, snap freezing in a dry/ice ethanol bath or liquid nitrogen. Freezing and storage of diluted enzyme is not recommended.



**Coomassie blue stained SDS-PAGE (4 – 12% acrylamide) of 4  $\mu$ g (right) of RBC PRMT6.** MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa



**Methyltransferase Activity of PRMT6.** Methylation determined as TCA-precipitable counts in a scintillation/filter plate assay. Reactions were 60 min., 30°C, with 1  $\mu$ M [<sup>3</sup>H]-SAM and 0.77 $\mu$ M protein substrates as indicated.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com