

## PRMT8 (GST/Strep) (Protein arginine N-methyltransferase 8; HRMT1L3, HRMT1L4)

**CATALOG NO.:** HMT-21-657

**LOT NO.:**

**DESCRIPTION:** Human recombinant PRMT8 (residues 61-394 (C-terminus); Genbank Accession # NM\_019854) expressed with N-terminal GST fusion protein and C-terminal Strep-tag, in Sf9 insect cells. MW = 67.09 kDa. PRMT8, 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 ( $\omega$ -monomethylation) and the transfer of a second methyl group to the same nitrogen, yielding asymmetric dimethylarginine (aDMA)<sup>1</sup>. Although highly homologous (>80% identity) to the major arginine methyltransferase PRMT1, distinctive characteristics of PRMT8 include its unique, regulatory 76 residue N-terminal domain and its brain-specific expression<sup>1</sup>, especially enhanced in the cortex<sup>2</sup>. Although transfected GFP-PRMT8 can be myristoylated and localized to the plasma membrane<sup>1</sup>, other evidence suggests that the endogenous, central nervous system PRMT8 is a nuclear protein<sup>3</sup>. Removal of the N-terminal 60 amino acids of PRMT8 has been shown to dramatically enhance PRMT8 catalytic activity<sup>4</sup>. RBC's PRMT8 is deleted for residues 1-60 and is highly active with the substrate GST-GAR.

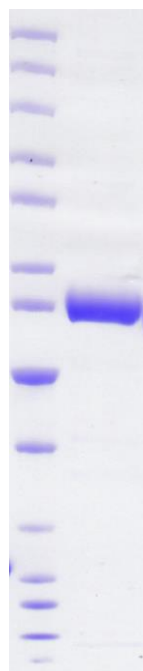
**PURITY:** >90% by SDS-PAGE.

**ASSAY CONDITIONS:** RBC's PRMT8 displays methyltransferase activity at enzyme concentrations of 15.6 nM and above, with GST-GAR (GST fused to fibrillarin residues 2-148) and [<sup>3</sup>H]-S-adenosyl-L-methionine as substrates. 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, 30°C, 60 min. with substrates as indicated above.

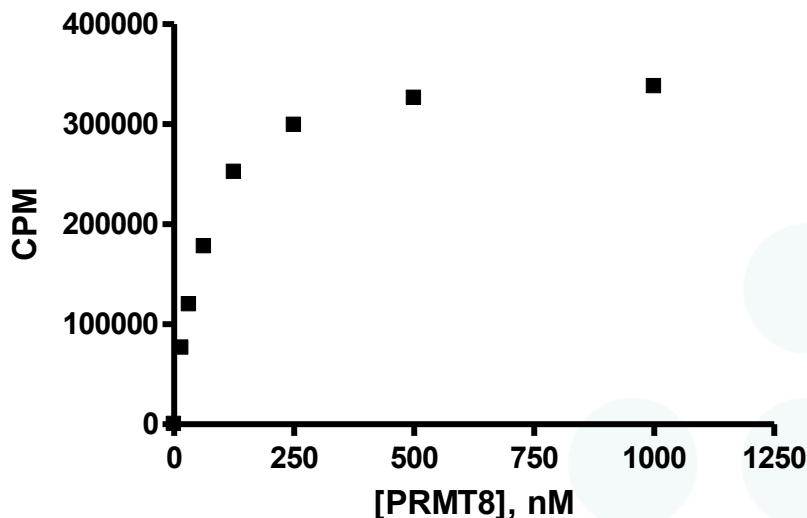
**SUPPLIED AS:** 2.13  $\mu$ g/ $\mu$ l in 50mM Tris-HCl pH 8, 500mM NaCl, 10% glycerol, 0.25mM TCEP, 2 mM DTT as determined by OD<sub>280</sub>

**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.

**REFERENCES:** 1) J. Lee *et al.* *J. Biol. Chem.* 2005 **280** 32890; 2) M.K. Weng *et al.* *PLOS One* 2012 **7** e36708; 3) A. Kousaka *et al.* *Neuroscience* 2009 **163** 1146; 4) J. Sayegh *et al.* *J. Biol. Chem.* 2007 **282** 36444



**Coomassie blue stained SDS-PAGE (4-12% acrylamide) of 4  $\mu$ g of RBC PRMT8.** MW markers (left) are, from top, 200, 150, 120, 100, 85, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**Methyltransferase Activity of PRMT8.** Methylation determined as TCA-precipitable counts in a scintillation/filter plate assay. Reactions were 20 $\mu$ L, 60 min., 30°C, with 1  $\mu$ M GST-GAR and 1 $\mu$ M [<sup>3</sup>H]-SAM as substrates.

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