

Recombinant PRMT1 protein

Catalog No: 31325

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 2 µg/µl

Source: Human

Buffer Contents: 10 µg recombinant PRMT1 protein supplied at a concentration of 2 µg/µl in 20 mM Tris-HCl, pH 7.3, 100 mM KCl, 0.2 mM EDTA and 20% glycerol.

Protein Details: PRMT1 (protein arginine N-methyltransferase 1) is a type 1 PRMT, capable of monomethylating and asymmetrically dimethylating arginine residues. PRMT1 specifically methylates arginine 3 on histone H4 (H4R3). Recombinant PRMT1 was expressed as the full-length (accession number NP_001527.3) human protein with an N-terminal poly-histidine tag.

Application Notes: Recombinant PRMT1 is suitable for use in methyltransferase assays. A recommended starting point for the methyltransferase assay is 50-200 ng of recombinant protein per reaction. The molecular weight of the protein is ~47 kDa.

Storage and Guarantee: This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.