

Recombinant PRMT3 protein

Catalog No: 31412, 31812

Expressed In: Baculovirus

Quantity: 20 µg

Concentration: 0.5 µg/µl

Source: Human

Buffer Contents: Full length recombinant PRMT3 protein is supplied at a concentration of 0.5 µg/µl in 25 mM HEPES pH 7.5, 120 mM NaCl, 5% glycerol, 0.2 mg/ml 3×FLAG peptide.

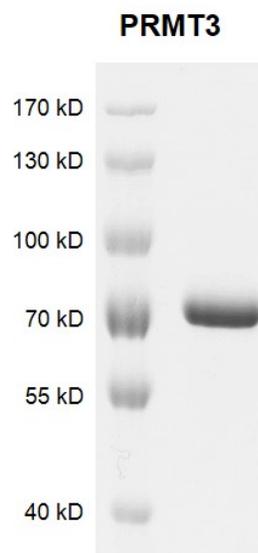
Background: PRMT3 (protein arginine N-methyltransferase 3) is a member of the protein arginine N-methyltransferase (PRMT) family that is capable of monomethylating and asymmetrically dimethylating arginine residues of histone H4 (H4R3). PRMT3 has been shown to interact with RPS2.

Protein Details: Recombinant PRMT3 (accession number NP_005779.1) was expressed in Sf9 cells and contains an N-terminal FLAG tag with an observed molecular weight of 61.5 kDa. The recombinant protein is >95% pure by SDS-PAGE.

Application Notes: Recombinant PRMT3 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. 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.



Recombinant PRMT3 protein gel. PRMT3 protein was run on a 8% SDS-PAGE gel and stained with Coomassie Blue.