

PRMT3 Polyclonal Antibody

(Catalog # A70419)

Background

This gene belongs to the protein arginine methyltransferase (PRMT) family. The encoded enzyme catalyzes the methylation of guanidino nitrogens of arginyl residues of proteins. The enzyme acts on 40S ribosomal protein S2 (rpS2), which is its major in-vivo substrate, and is involved in the proper maturation of the 80S ribosome. Alternative splicing results in multiple transcript variants.

Description

PRMT3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

O60678

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-220 of human PRMT3 (NP_005779.1).

Storage

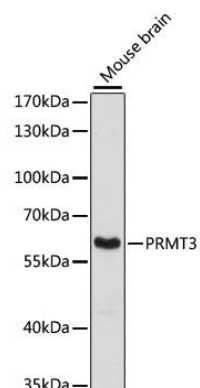
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

PRMT3; HRMT1L3; protein arginine methyltransferase 3

Application

WB; Recommended dilution: WB 1:1000 - 1:2000



Western blot analysis of extracts of mouse brain, using PRMT3 Polyclonal Antibody at 1:3000 dilution.
Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 90s.