
PRMT1 Polyclonal Antibody

(Catalog # A-3001)

Background

PRMTs (Protein Arginine N-Methyltransferases) can monomethylate or dimethylate arginine residues by transferring methyl groups from S-adenoxylmethionine to the guanidine nitrogen of arginine. PRMT1, also known as Protein Arginine N-Methyltransferase 1, has been shown to play a role in transcriptional regulation, specifically methylating Histone H4R3 and acting alongside p300/CBP. It is also able to methylate non-histone proteins. PRMT1 is implicated in other activities such as signal transduction, mRNA processing, and p53-mediated transcription.

Description

PRMT1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse, Rat, Human

Isotype

IgG

Uniprot ID

Q99873

Purification

Affinity Purified

Immunogen

Recombinant Protein of Human PRMT1

Storage

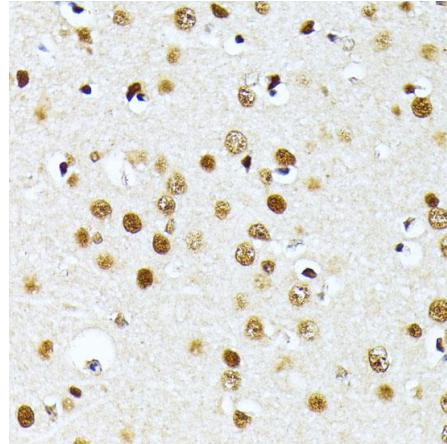
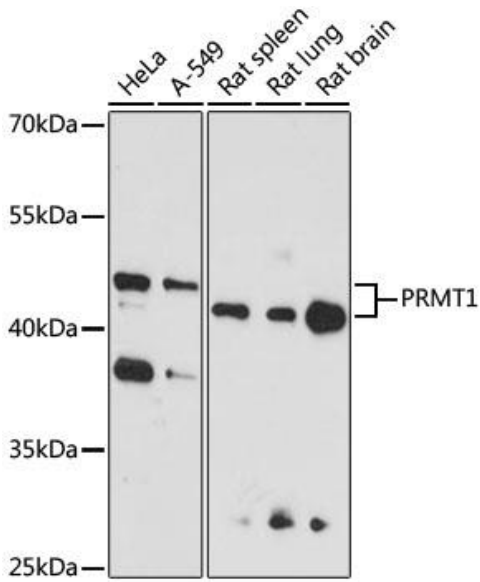
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

ANM 1 antibody, ANM1_HUMAN antibody, complete cds antibody, HCP 1 antibody, Heterogeneous nuclear ribonucleoprotein methyltransferase 1 like 2 antibody, Histone-arginine N-methyltransferase PRMT1 antibody, HMT 2 antibody, HMT1 (hnRNP methyltransferase *S. cerevisiae*) like 2 antibody, HMT1 hnRNP methyltransferase antibody, HMT2 antibody, HRMT1 L2 antibody, Human mRNA for suppressor for yeast mutant complete cds antibody, Interferon receptor 1 bound protein 4 antibody, IR1 B4 antibody, Mrmt 1 antibody, PRMT 1 antibody, Protein arginine methyltransferase 1 antibody, Protein arginine N-methyltransferase 1 (EC 2.1.1.-) (Interferon receptor 1-bound protein 4) antibody, R1B4 antibody

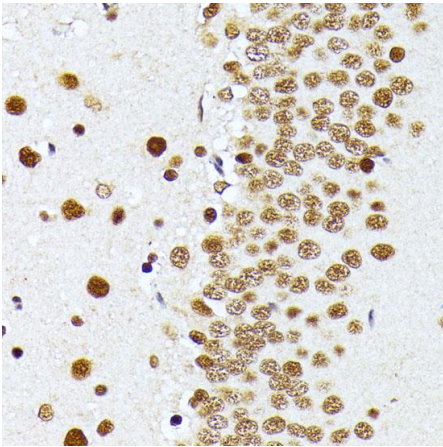
Application

WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000; IHC 1:50 - 1:200; IF 1:50 - 1:200

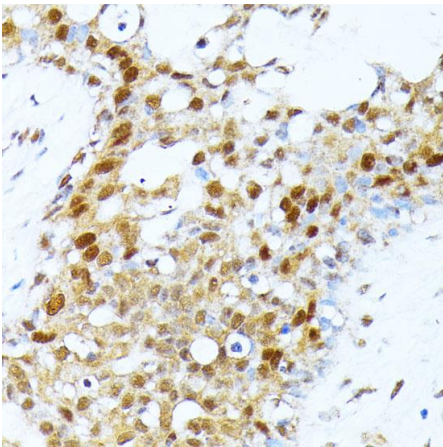


Immunohistochemistry of paraffin-embedded mouse brain using PRMT1 antibody at dilution of 1:100 (40x lens).

Western blot analysis of extracts of various cell lines, using PRMT1 antibody at 1:1000 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.



Immunohistochemistry of paraffin-embedded rat brain using PRMT1 antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human breast cancer using PRMT1 antibody at dilution of 1:100 (40x lens).