
PRMT4 Polyclonal Antibody

(Catalog # A-3004)

Background

PRMT4 (Protein Arginine N-Methyltransferase 4), also known as CARM1 (coactivator-associated arginine methyltransferase 1), is a part of the PRMT family that can monomethylate or dimethylate arginine residues by transferring methyl groups from S-adenoxylmethionine to the guanidine nitrogen of arginine. PRMT4 has been shown to catalyze methylation at histone H3 Arg-17 (H3R17) and Histone 3 Arg-26 (H3R26).

Description

PRMT4 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

PBS with 0.05% proclin300, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q86X55

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 486-603 of human PRMT4/CARM1 (NP_954592.1)

Storage

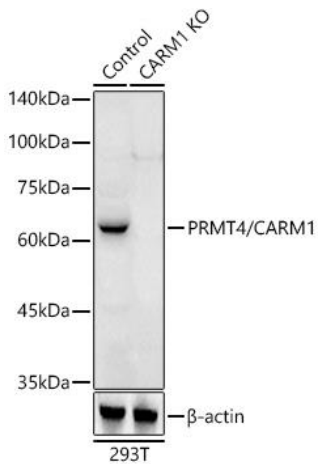
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze.

Alternative Names

carm1 antibody, CARM1_HUMAN antibody, Coactivator associated arginine methyltransferase 1 antibody, Histone arginine methyltransferase CARM 1 antibody, PRMT 4 antibody, Protein arginine N methyltransferase 4 antibody

Application

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

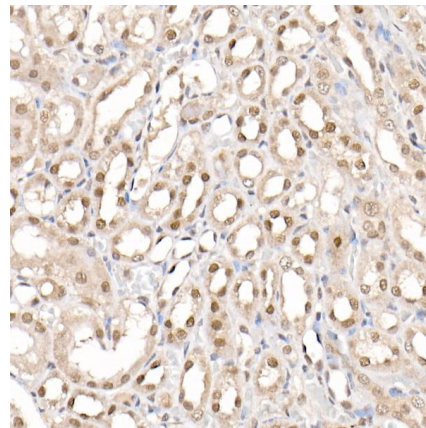


Western blot analysis of extracts from normal (control) and PRMT4/CARM1 knockout (KO) 293T cells, using PRMT4 Polyclonal Antibody at 1:500 dilution.

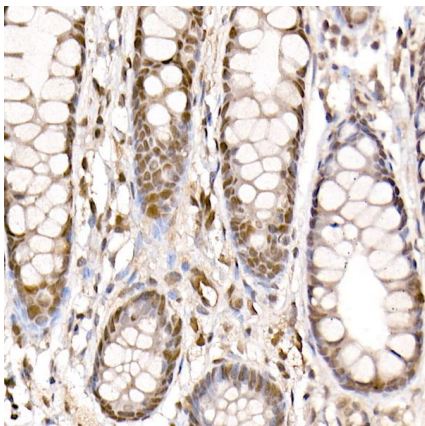
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

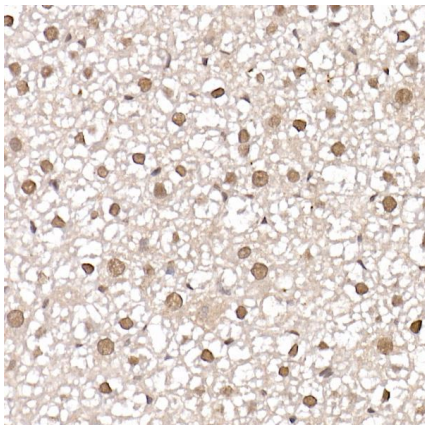
Blocking buffer: 3% nonfat dry milk in TBST.



Immunohistochemistry of paraffin-embedded rat kidney using PRMT4 Polyclonal Antibody at dilution of 1:25 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry of paraffin-embedded human colon using PRMT4 Polyclonal Antibody at dilution of 1:25 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry of paraffin-embedded mouse liver using PRMT4 Polyclonal Antibody at dilution of 1:25 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.