
PRMT6 Polyclonal Antibody

(Catalog # A70420)

Background

The protein encoded by this gene belongs to the arginine N-methyltransferase family, which catalyze the sequential transfer of methyl group from S-adenosyl-L-methionine to the side chain nitrogens of arginine residues within proteins, to form methylated arginine derivatives and S-adenosyl-L-homocysteine. This protein can catalyze both, the formation of omega-N monomethylarginine and asymmetrical dimethylarginine, with a strong preference for the latter. It specifically mediates the asymmetric dimethylation of Arg2 of histone H3, and the methylated form represents a specific tag for epigenetic transcriptional repression. This protein also forms a complex with, and methylates DNA polymerase beta, resulting in stimulation of polymerase activity by enhancing DNA binding and processivity.

Description

PRMT6 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thiomersal, 50% glycerol, pH7.3.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q96LA8

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-375 of human PRMT6 (NP_060607.2).

Storage

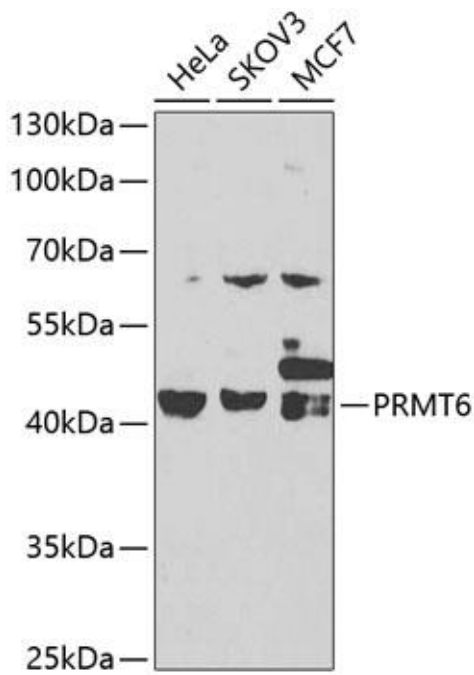
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

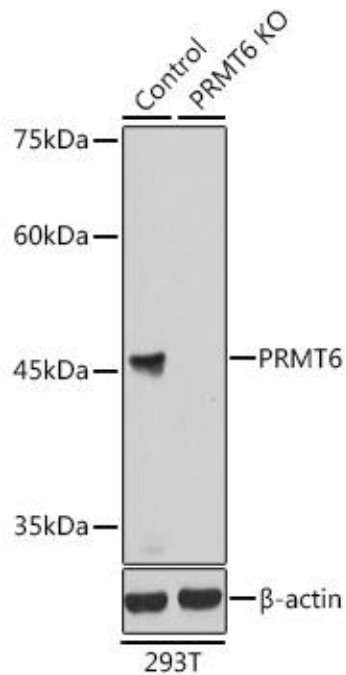
PRMT6; HRMT1L6; protein arginine methyltransferase 6

Application

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



Western blot analysis of extracts of various cell lines, using PPRMT6 Polyclonal Antibody at 1:1000 dilution.
 Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25µg per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.



Western blot analysis of extracts from normal (control) and PRMT6 Rabbit pAb knockout (KO) 293T cells, using PPRMT6 Polyclonal Antibody at 1:1000 dilution.
 Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25µg per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.