

## PRMT6 Polyclonal Antibody

(Catalog # A-3006)

---

### Background

PRMT6, or Protein arginine N-methyltransferase 6, is part of the PRMT family that catalyzes the methylation of arginine residues. PRMT6 specifically coregulates gene expression by acting as an activator and repressor via dimethylation of Histone H3R2me2a. PRMT6 methylates several histones at various residues, particularly Histone H2AR3me, Histone H3K4me2, Histone H3K4me3, as well as Histone H4R3me. By transcriptionally repressing tumor suppressor genes, PRMT6 has been shown to regulate cell proliferation and senescence and, via interactions with Rev, Nucleocapsid protein p7, and Tat, PRMT6 might also be implicated in innate immunity against HIV-1.

### Description

PRMT6 Polyclonal Antibody. Unconjugated. Raised in: Rabbit

### Formulation

Liquid. 0.03% Proclin 300. 50% Glycerol, 0.01M PBS, PH 7.4

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q96LA8

### Purification

Protein G purified

### Immunogen

Recombinant Human Protein arginine N-methyltransferase 6 protein (119-375AA)

### Storage

Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

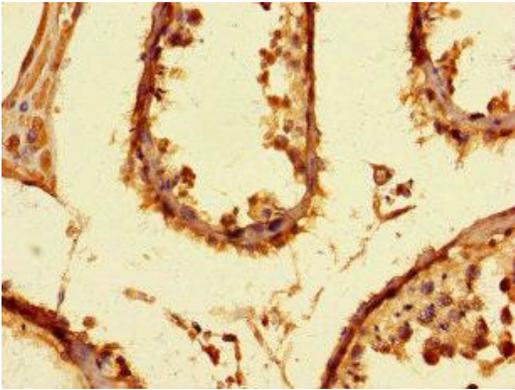
### Alternative Names

Protein arginine N-methyltransferase 6 (EC 2.1.1.319) (Heterogeneous nuclear ribonucleoprotein methyltransferase-like protein 6) (Histone-arginine N-methyltransferase PRMT6), PRMT6, HRMT1L6

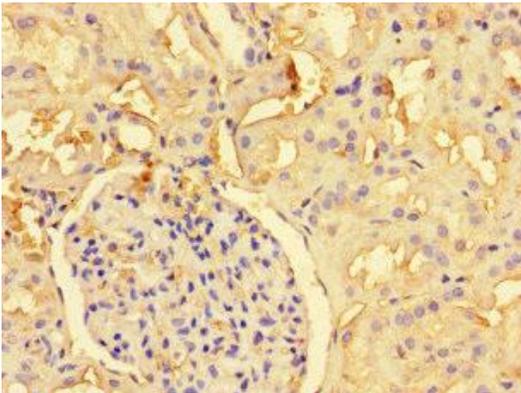
### Application

ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200

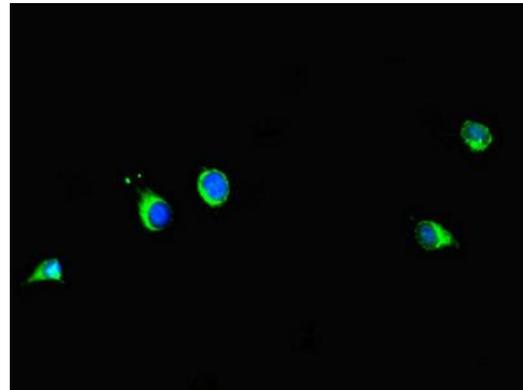
---



Immunohistochemistry of paraffin-embedded human testis tissue using PRMT6 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human kidney tissue using PRMT6 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HeLa cells using PRMT6 Polyclonal Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)