

PRMT5 antibody (pAb)

Catalog Nos: 61001, 61002

RRID: AB_2615010

Isotype: IgG

Application(s): ChIP-Seq, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.65 µg/µl

Molecular Weight: 75 kDa

Background: PRMT5 (Protein Arginine Methyltransferase 5, also known as JBP1, Janus Binding Protein 1) is a type II arginine methyltransferase, a protein that transfers methyl groups to the arginine residues of substrate proteins. Arginine methylation is a common post-translational modification of histones and other cellular proteins. PRMT5 symmetrically dimethylates histone H2A Arg3, H3 Arg8 and H4 Arg3. It also methylates PIWI proteins which regulate small non-coding RNAs and snRNPs, regulating snRNP core particle assembly. PRMT5 functions as a co-repressor of transcription and is recruited by a number of DNA binding proteins and transcriptional repressor complexes. It is present both in the nucleus and the cytoplasm. It also interacts with COPR5, JAK2, PRDM1 and RPS10.

Immunogen: This PRMT5 antibody was raised against a peptide corresponding to amino acids within the N-terminal region of human PRMT5.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

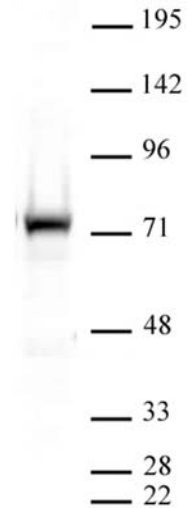
WB: 1:500 - 1:2,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

References: PRMT5 antibody has been cited in one or more publications. To see an up-to-date listing of papers that describe its use, please go to www.activemotif.com/catalog/details/61001.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Western blot of PRMT5 antibody.

Cytoplasmic extract of Jurkat cells (30 µg per lane) probed with PRMT5 antibody (1:1,000).