

## PRMT2 antibody (pAb)

**Catalog Nos:** 61491, 61492

**RRID:** AB\_2793656

**Isotype:** IgG

**Application(s):** ICC, IF, WB

**Reactivity:** Human

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 55 kDa

**Background:** PRMT2 (Protein Arginine Methyltransferase 2) is a member of the protein arginine N-methyltransferase (PRMT) family that is capable of monomethylating and asymmetrically dimethylating arginine residues. PRMT2 is an arginine methyltransferase that methylates the guanidino nitrogens of arginyl residues in proteins such as STAT3, FBL, histone H4. Acts as a coactivator (with NCOA2) of the androgen receptor (AR)-mediated transactivation. Acts as a coactivator (with estrogen) of estrogen receptor (ER)-mediated transactivation. Enhances PGR, PPAR $\gamma$ , RARA-mediated transactivation. May inhibit NF-kappa-B transcription and promote apoptosis. Represses E2F1 transcriptional activity (in a RB1-dependent manner). May be involved in growth regulation.

**Immunogen:** This antibody was raised against a peptide within the N-terminal region of human PRMT2.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

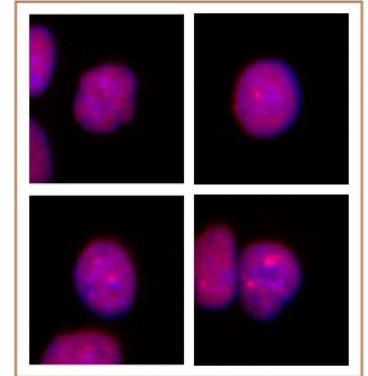
Applications Validated by Active Motif:

ICC/IF: 1:2,000 dilution

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

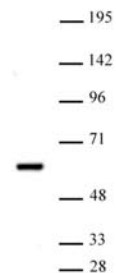
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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



### PRMT2 antibody (pAb) tested by Immunofluorescence.

Formaldehyde fixed HeLa cells stained with PRMT2 antibody at a 1:2,000 dilution.



### PRMT2 antibody (pAb) tested by Western blot.

Nuclear extract of U373 MG cells (20  $\mu$ g) probed with PRMT2 antibody (pAb) at a dilution of 1:1,000.