

PRMT7 polyclonal antibody - Classic

Cat. No. C15410247

Type: Polyclonal

Source: Rabbit

Lot #: 40674

Size: 25 µl/100 µl

Concentration: 1 µg/µl

Specificity: Human, mouse, rat.

Purity: Purified by antigen-affinity chromatography.

Storage: Store at -20°C; for long storage, store at -80°C.
Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Applications

	Suggested dilution*	Results
ChIP assay	5 µg/ChIP	Figure 1
ICC/IF	1:100-1:1,000	Figure 3
IHC	1:100-1:1,000	Figure 5, 6, 7
Immunoprecipitation	1:100-1:500	Figure 2
Western blot	1:500-1:3,000	Figure 4

*Optimal dilutions/concentrations should be determined by the researcher.

Results

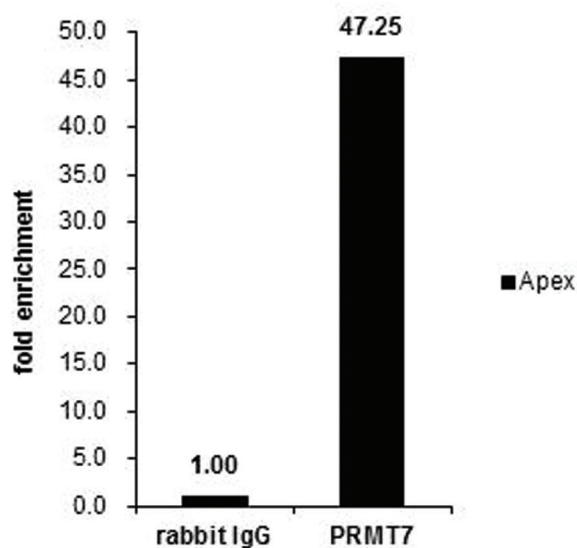


Figure 1. ChIP

ChIP was performed with NIH3T3 chromatin extract and 5 µg of either control rabbit IgG or PRMT7 antibody. The precipitated DNA was detected by PCR with primer set targeting to Apex2.

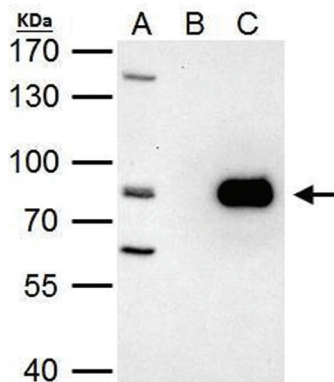


Figure 2. Immunoprecipitation

PRMT7 antibody immunoprecipitates PRMT7 protein in IP experiments. IP Sample: HeLa whole cell extract A. 40 µg HeLa whole cell extract B. Control with 2 µg of preimmune rabbit IgG C. Immunoprecipitation of PRMT7 protein by 2 µg of PRMT7 antibody (Cat. No. C15410247) The immunoprecipitated PRMT7 protein was detected by PRMT7 antibody (C15410247) diluted 1:1,000.

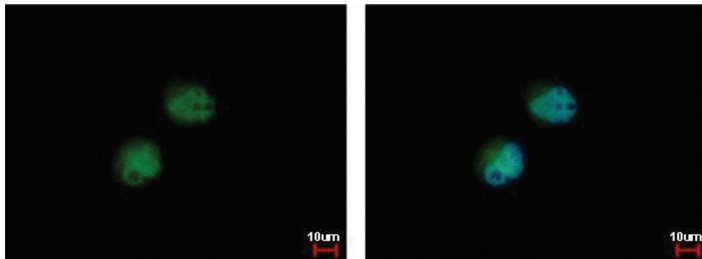


Figure 3. Immunofluorescent analysis

PRMT7 antibody detects PRMT7 protein in the cytoplasm and nucleus by immunofluorescent analysis.

Sample: HeLa cells were fixed in 4% paraformaldehyde at RT for 15 min.

Green: PRMT7 protein stained by PRMT7 antibody (Cat. No. C15410247) diluted 1:500.

Blue: Hoechst 33342 staining.

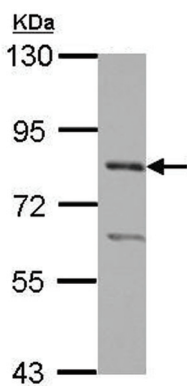


Figure 4. Western blot

Sample: 30 µg of PC-3 whole cell lysate

7.5% SDS PAGE

PMRT7 antibody diluted 1:2,000

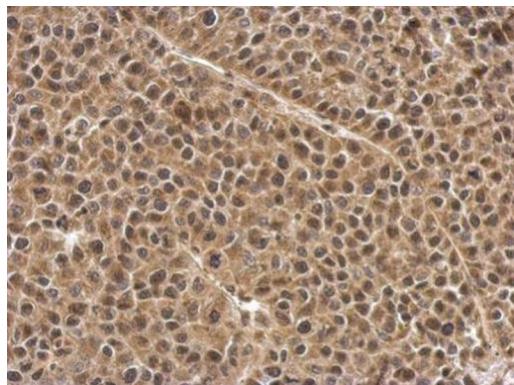


Figure 5. Immunohistochemical analysis

Immunohistochemical analysis of paraffin-embedded Hela xenograft, using PRMT7 (C15410247) antibody at a 1:500 dilution.

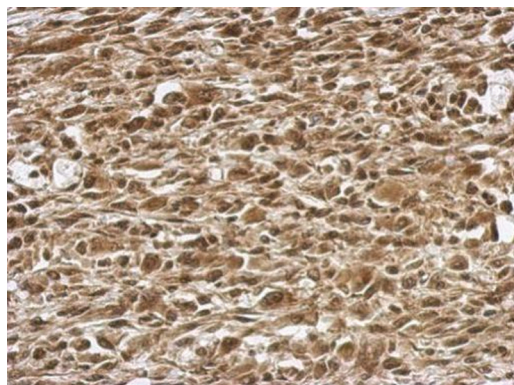


Figure 6. Immunohistochemical analysis

Immunohistochemical analysis of paraffin-embedded C2C12 xenograft, using PRMT7 (C15410247) antibody at a 1:500 dilution.

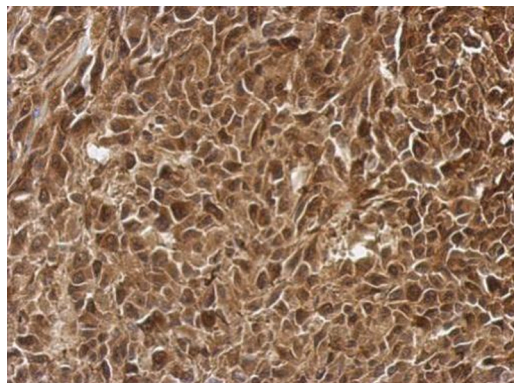


Figure 7. Immunohistochemical analysis

Immunohistochemical analysis of paraffin-embedded RT2 xenograft, using PRMT7 (C15410247) antibody at a 1:500 dilution.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: September 12, 2014