

DNMT1 polyclonal antibody - Classic

Cat. No. C15410227

Type: Polyclonal

Source: Rabbit

Lot #: 40135

Size: 25 µl/100 µl

Concentration: 0,74 µg/µl

Specificity: Human.

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 6
ICC/IF	1:100-1:1,000	Figure 2
IHC	1:100-1:1,000	Figure 1
Immunoprecipitation	1:100-1:500	Figure 5
Western blot	1:500-1:3,000	Figure 3, 4

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

Results

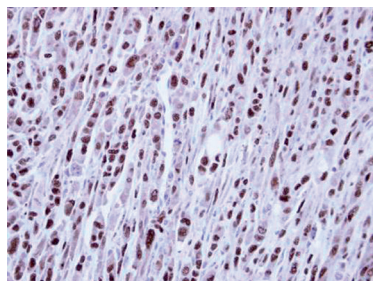


Figure 1. IHC

Immunohistochemical analysis of paraffin-embedded PC3 Xenograft, using DNMT1 (C15410227) antibody at a 1:100 dilution.

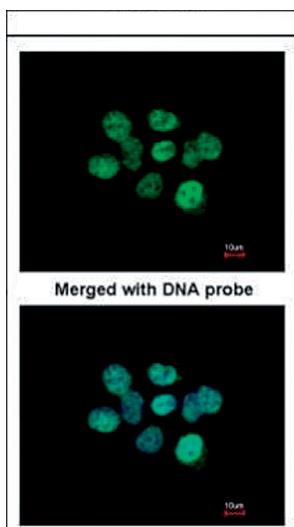


Figure 2. IF

Immunofluorescence analysis of paraformaldehyde-fixed A431, using DNMT1 (C15410227) antibody at a 1:200 dilution.

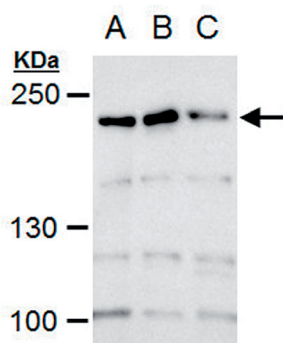


Figure 3. Western blot

DNMT1 antibody detects DNMT1 protein by western blot analysis.

A. 30 μ g 293T whole cell extract
B. 30 μ g A431 whole cell extract
C. 30 μ g HeLa whole cell extract
D. 30 μ g HepG2 whole cell extract
5 % SDS-PAGE
DNMT1 antibody (C15410227) dilution: 1:500

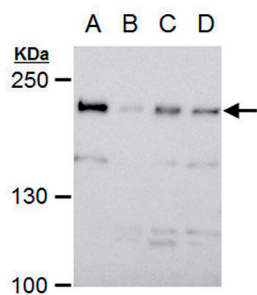


Figure 4. Western blot

DNMT1 antibody detects DNMT1 protein by western blot analysis.

A. 30 μ g Jurkat whole cell extract
B. 30 μ g Raji whole cell extract
C. 30 μ g NCI-H929 whole cell extract
5 % SDS-PAGE
DNMT1 antibody (C15410227) dilution: 1:500

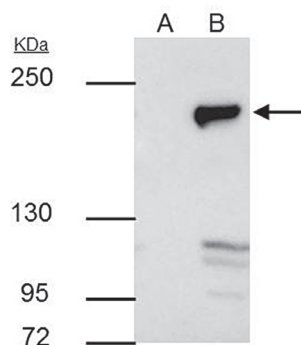


Figure 5. IP

DNMT1 antibody immunoprecipitates DNMT1 protein in IP experiments.

IP samples: HeLa nuclear extract

A. Control with 4 µg of preimmune Rabbit IgG

B. Immunoprecipitation of DNMT1 protein by 4 µg DNMT1 antibody (C15410227)

5 % SDS-PAGE

The immunoprecipitated DNMT1 protein was detected by DNMT1 antibody (C15410227) diluted 1:1,000.

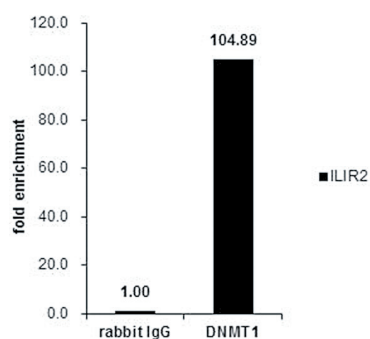


Figure 6. ChIP

ChIP was performed with HCT116 chromatin extract and 5 µg of either control rabbit IgG or DNMT1 antibody. The precipitated DNA was detected by PCR with primer set targeting to ILIR2. The results are expressed as fold enrichment over the IgG negative control.

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