

## 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.

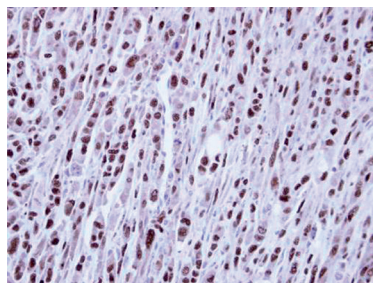
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

### 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

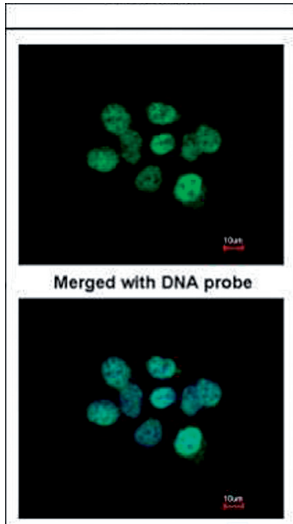
\*Optimal dilutions/concentrations should be determined by the researcher.

### Results



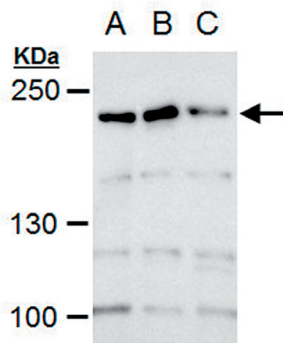
**Figure 1. IHC**

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



**Figure 2. IF**

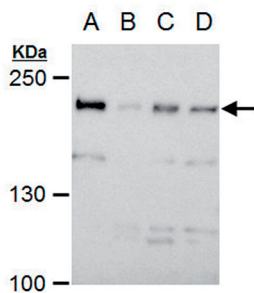
Immunofluorescence analysis of paraformaldehyde-fixed A431, using DNMT1 (C15410227) antibody at a 1:200 dilution.



**Figure 3. Western blot**

DNMT1 antibody detects DNMT1 protein by western blot analysis.

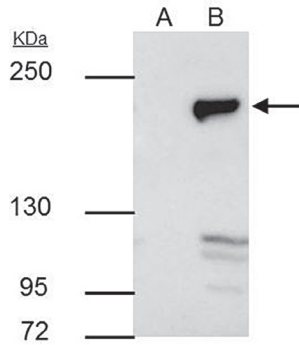
- A. 30 µg 293T whole cell extract
  - B. 30 µg A431 whole cell extract
  - C. 30 µg HeLa whole cell extract
  - D. 30 µg HepG2 whole cell extract
- 5 % SDS-PAGE  
DNMT1 antibody (C15410227) dilution: 1:500



**Figure 4. Western blot**

DNMT1 antibody detects DNMT1 protein by western blot analysis.

- A. 30 µg Jurkat whole cell extract
  - B. 30 µg Raji whole cell extract
  - C. 30 µg NCI-H929 whole cell extract
  - D. 30 µg HepG2 whole cell extract
- 5 % SDS-PAGE  
DNMT1 antibody (C15410227) dilution: 1:500



**Figure 5. IP**

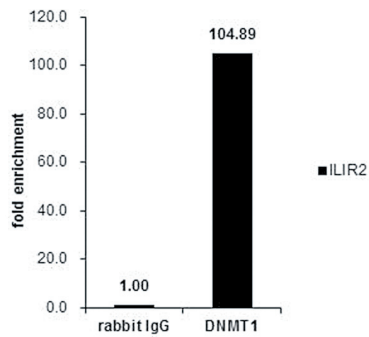
DNMT1 antibody immunoprecipitates DNMT1 protein in IP experiments.

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