

SATB1 polyclonal antibody - Classic

Cat. No. C15410248

Type: Polyclonal

Source: Rabbit

Lot #: 40499

Size: 25 µl/100 µl

Concentration: 0.81 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 1
IHC	1:100-1:1,000	Figure 4
Immunoprecipitation	1:100-1:500	Figure 2
Western blot	1:500-1:3,000	Figure 3, 5

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

Results

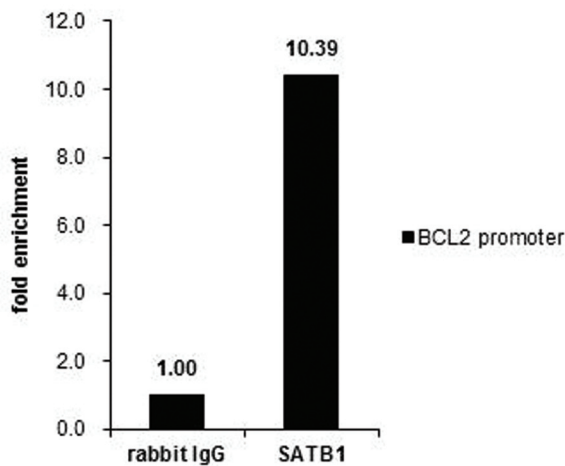


Figure 1. ChIP

ChIP was performed with Jurkat chromatin extract and 5 µg of either control rabbit IgG or SATB1 antibody. The precipitated DNA was detected by PCR with primer set targeting to BCL2 promoter.

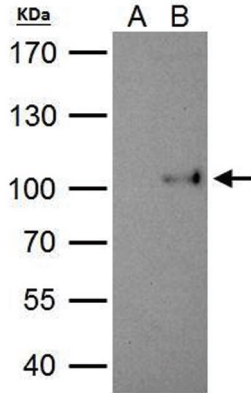


Figure 2. Immunoprecipitation

SATB1 antibody immunoprecipitates SATB1 protein in IP experiments. IP Sample: 293T whole cell extract.

A. Control with 2 μ g of preimmune rabbit IgG (Cat. No. C15410206).

B. Immunoprecipitation of SATB1 protein by 2 μ g of SATB1 antibody
7.5% SDS-PAGE

The immunoprecipitated SATB1 protein was detected by SATB1 antibody diluted 1:1,000.

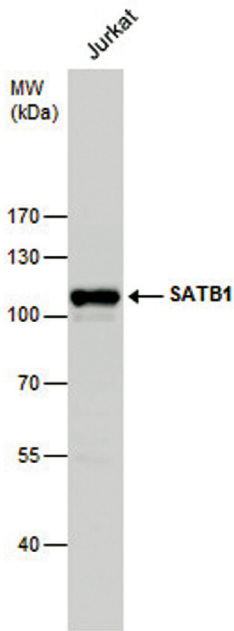


Figure 3. Western Blot

SATB1 antibody detects SATB1 protein by western blot analysis. Jurkat whole cell extracts (30 μ g) were separated by 7.5 % SDS-PAGE, and blotted with SATB1 antibody (C15410248) diluted 1:500

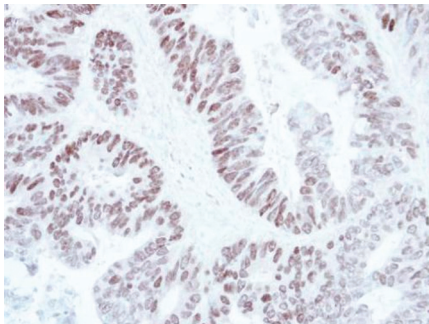


Figure 4. Immunohistochemistry

Immunohistochemical analysis of paraffin-embedded Colon, using the SATB1 antibody (C15410248) at a 1:250 dilution.

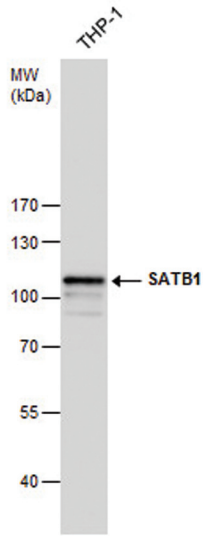


Figure 5. Western Blot

SATB1 antibody detects SATB1 protein by western blot analysis. THP-1 whole cell extracts (30 µg) were separated by 7.5 % SDS-PAGE, and blotted with SATB1 antibody (C15410248) diluted 1:500

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