

ZEB1 polyclonal antibody - Classic

Cat. No. C15410259

Type: Polyclonal

Source: Rabbit

Lot #: 42690

Size: 25 µl /100 µl

Concentration: 0.1 µg / µl

Specificity: Human.

Purity: Purified by peptide 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 4
Immunoprecipitation	1:100-1:500	Figure 2
Western blot	1:500-1:3,000	Figure 3

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

Results

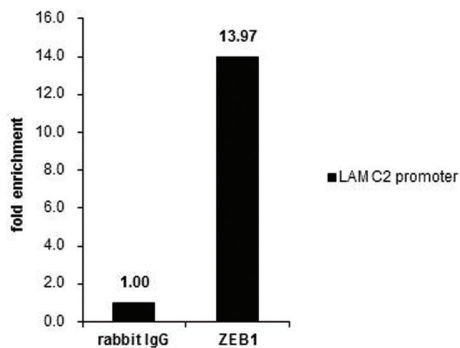


Figure 1. ZEB1 ChIP results

ChIP was performed with HeLa chromatin extract and 5 µg of either control rabbit IgG or ZEB1 antibody. The precipitated DNA was detected by PCR with primer set targeting to LAMC2 promoter.

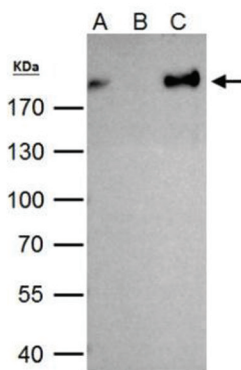


Figure 2. IP results

ZEB1 antibody immunoprecipitates ZEB1 protein in IP experiments. IP Sample: 293T whole cell extract A. 40 µg 293T whole cell extract B. Control with 2 µg of preimmune rabbit IgG C. Immunoprecipitation of ZEB1 protein by 2 µg of ZEB1 antibody (Cat. No. C15410259). The immunoprecipitated ZEB1 protein was detected by western blot with the ZEB1 antibody diluted 1:1,000.

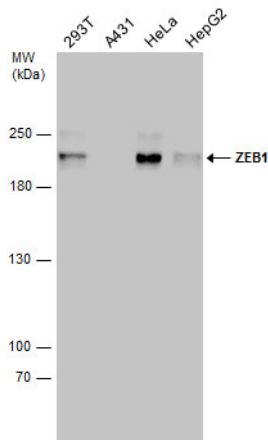


Figure 3. Western blot

Whole cell extracts (30 μ g) from 293T, A431, HeLa and HepG2 cells were analyzed by WB using the Diagenode ZEB1 antibody (C15410259) diluted 1:1000. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

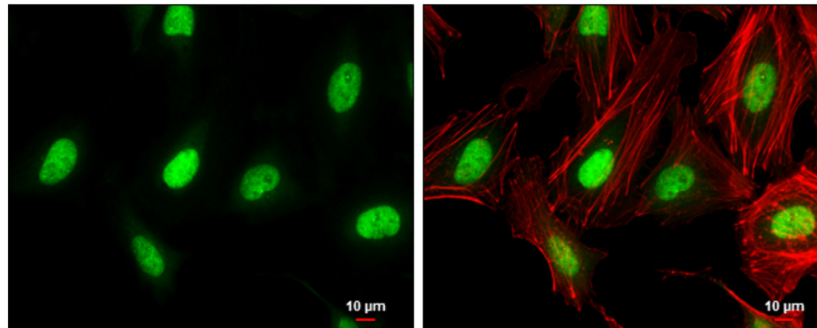


Figure 4. IF

Immunofluorescence analysis of paraformaldehyde-fixed HeLa cells using the Diagenode ZEB1 antibody (Cat. No. C15410259) at a 1:500 dilution (green). The cytoskeleton was stained with phalloidin (red).

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: March 27, 2017