

CTNNB1 polyclonal antibody - Classic

Other names: Catenin Beta, MRD19

Cat. No. C15410222

Type: Polyclonal ChIP-grade

Source: Rabbit

Lot #: 42235

Size: 25 µl/100 µl

Concentration: 0.1 µg/µl

Specificity: Human, mouse, rat: positive

Other species: not tested

Purity: Affinity purified polyclonal antibody in PBS containing 1% BSA, 20% Glycerol and 0.025% ProClin 300.

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.

Description: Polyclonal antibody raised in rabbit against CTNNB1 (Catenin Beta 1), using a recombinant protein.

Applications

	Suggested dilution/amount	Results
ChIP*	5 µg per ChIP	Fig 1
Western blotting	1:1,000 - 1:3,000	Fig 2, 3
Immunoprecipitation	5 µg per IP	Fig 4
Immunofluorescence	1:500	Fig 5
Immunohistochemistry	1:500	Fig 6

* Please note that the optimal antibody amount per IP should be determined by the end-user. We recommend testing 1-5 µg per IP.

Target description

CTNNB1 (UniProt/Swiss-Prot entry P35222) is a component of the Wnt signalling pathway. In the presence of Wnt, it accumulates in the nucleus where it acts as a coactivator for transcription factors of the TCF/LEF family, leading to the activation of Wnt responsive genes. CTNNB1 is also part of a protein complex that constitutes adherens junctions which are necessary for the creation and maintenance of epithelial cell layers by regulating cell growth and cell adhesion. Mutations in the CTNNB1 gene are a cause of colorectal cancer, pilomatixoma, medulloblastoma and ovarian cancer.

Results

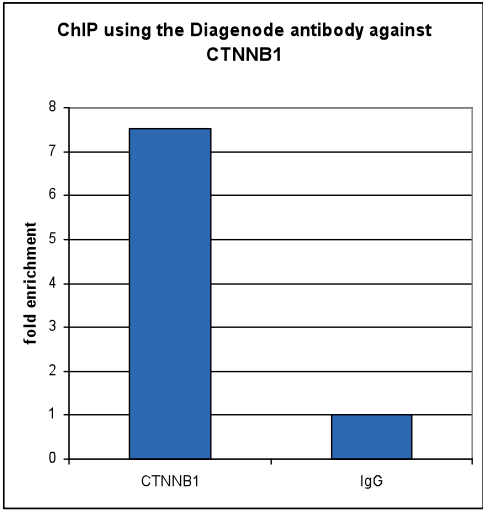


Figure 1. ChIP results obtained with the Diagenode antibody directed against CTNNB1

ChIP assays were performed using HCT116 cells and 5 µg of the Diagenode antibody against CTNNB1 (Cat. No. C15410222). An equal amount of IgG was used as a negative IP control. QPCR was performed with primers specific for the c-Myc promoter. Figure 1 shows the fold enrichment over the IgG negative control.

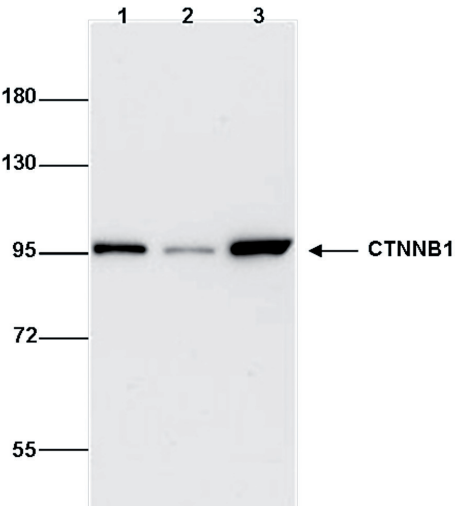


Figure 2. Western blot analysis using the Diagenode antibody directed against CTNNB1

Whole cell extracts (30 µg) from A549 (lane1), H1299 (lane 2) and HCT116 (lane 3) cells were analysed by Western blot using the Diagenode antibody against CTNNB1 (Cat. No. C15410222) diluted 1:3,000. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

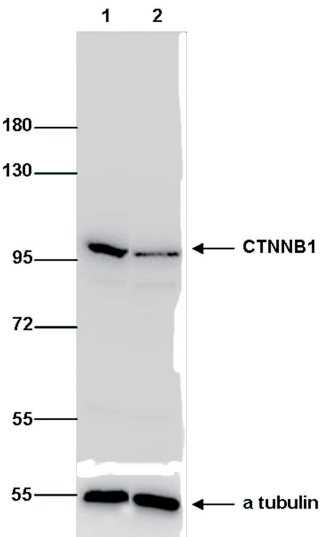


Figure 3. Western blot analysis using the Diagenode antibody directed against CTNNB1

Whole cell extracts (30 µg) from HeLa cells transfected with CTNNB1 shRNA (lane 2) and from an untransfected control (lane 1) were analysed by Western blot using the Diagenode antibody against CTNNB1 (Cat. No. C15410222) diluted 1:1,000. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left. Alpha tubulin was used as a loading control (lower panel).

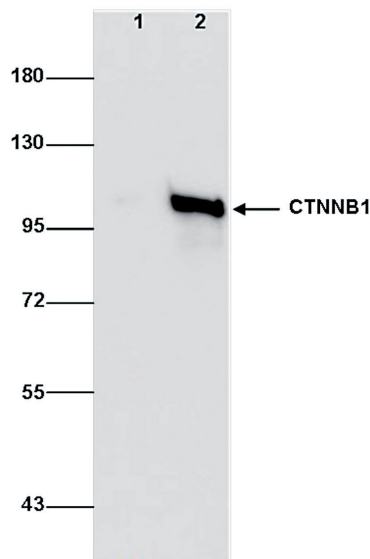


Figure 4. Immunoprecipitation using the Diagenode antibody directed against CTNNB1

Immunoprecipitation was performed on whole cell extracts from HeLa cells using 5 µg of the Diagenode antibody against CTNNB1 [Cat. No. C15410222, lane 2]. An equal amount of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated CTNNB1 protein was detected by western blot with the CTNNB1 antibody diluted 1:1,000.

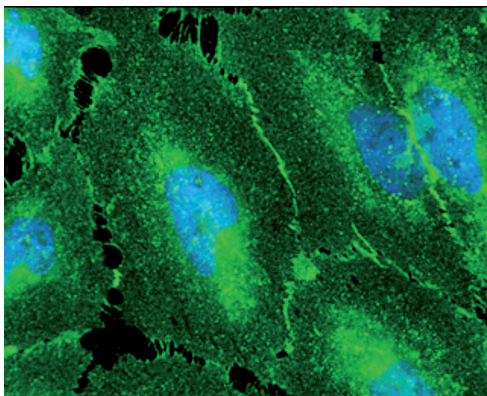


Figure 5. Immunofluorescence with the Diagenode antibody directed against CTNNB1

HeLa cells were fixed with 4% formaldehyde for 15' at room temperature and stained with the Diagenode antibody against CTNNB1 [Cat. C15410222] diluted 1:500 (green). The cell nuclei were stained with Hoechst 33342 (blue).

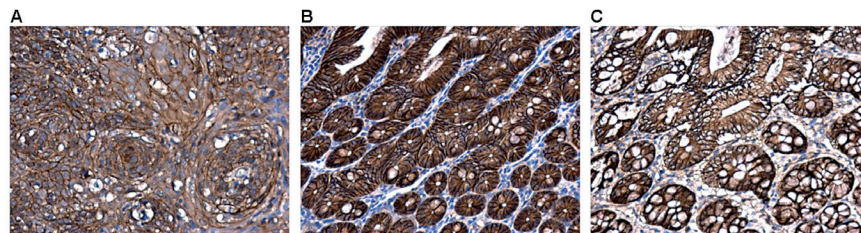


Figure 6. Immunohistochemistry using the Diagenode antibody directed against CTNNB1

Formalin fixed paraffin embedded human cervix (figure A), mouse intestine (figure B) and rat colon (figure C) were stained with the Diagenode antibody against CTNNB1 [Cat. No. C15410222] diluted 1:500, followed by a peroxidase labelled goat anti-rabbit secondary antibody.

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: October 18, 2016