

Brn2 polyclonal antibody - Classic

Cat. No. C15410223

Type: Polyclonal

Source: Rabbit

Lot #: 40212

Size: 25 µg/100 µg

Concentration: 1 µg/µl

Specificity: Human, mouse.

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
ICC/IF	1:100-1:1,000	
IHC	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

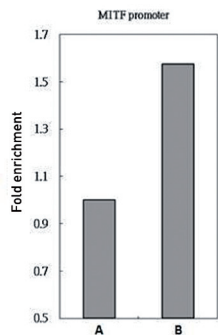


Figure 1. ChIP

Brn2 antibody immunoprecipitates Brn2 protein-DNA in ChIP experiments. ChIP Sample: 293T chromatin extract A. 5 µg preimmune rabbit IgG B. 5 µg of Brn2 antibody (C15410223) The precipitated DNA was detected by PCR with primer set targeting to MITF promoter.

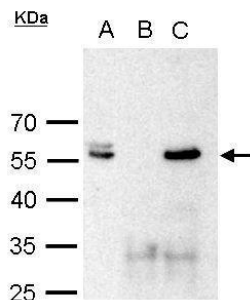


Figure 2. IP

Brn2 antibody immunoprecipitates Brn2 protein in IP experiments. IP Sample: 1000 µg 293T whole cell extract A. 50 µg 293T whole cell extract B. Control with 2 µg of preimmune rabbit IgG C. Immunoprecipitation of Brn2 protein by 2 µg of Brn2 antibody (C15410223) The immunoprecipitated Brn2 protein was detected by 10% SDS-PAGE using the Brn2 antibody (C15410223) diluted 1:1,000.

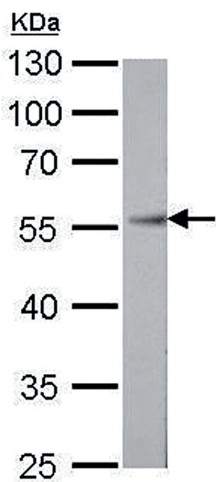


Figure 3. Western blot

Brn2 antibody detects Brn2 protein by Western blot analysis.
30 µg A375 whole cell extract
10 % SDS-PAGE
Brn2 antibody [C15410223] dilution: 1:1,000

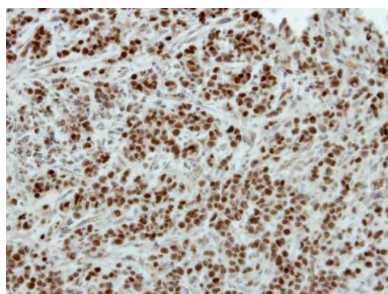


Figure 4. IHC

Immunohistochemical analysis of paraffin-embedded TOV-21G xenograft, using Brn2 [C15410223] antibody at a 1:500 dilution.

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 9, 2014