

DP1 polyclonal antibody - Classic

Cat. No. C15410228

Type: Polyclonal

Source: Rabbit

Lot #: 40422

Size: 25 µl/100 µl

Concentration: 1 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
Immunoprecipitation	1:100-1:500	Figure 1
Western blot	1:5,000-1:20,000	Figure 2

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

Results

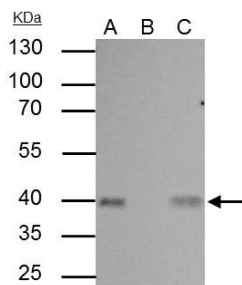


Figure 1. DP1 antibody IP results

DP1 antibody immunoprecipitates DP1 protein in IP experiments. IP Sample: 293T whole cell extract A : 40 µg whole cell extract of DP1 protein expressing 293T cells B : Control with 2.5 µg of pre-immune rabbit IgG (Cat. No. C15410206) C : Immunoprecipitation of DP1 by 2.5 µg of DP1 antibody (Cat. No. C15410228) 10% SDS-PAGE The immunoprecipitated DP1 protein was detected by DP1 antibody (Cat. No. C15410228) diluted 1:1,000.

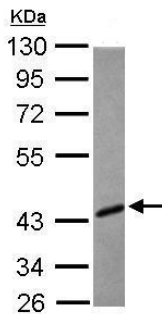


Figure 2. Western blot

Sample: 30 µg of whole cell lysate (raji cell)
10% SDS PAGE
C15410228 diluted 1:10,000

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 10, 2014