

UHRF1 polyclonal antibody - Classic

Cat. No. C15410258

Type: Polyclonal

Source: Rabbit

Lot #: 43922

Size: 25 µl/100 µl

Concentration: 0.97 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 5
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

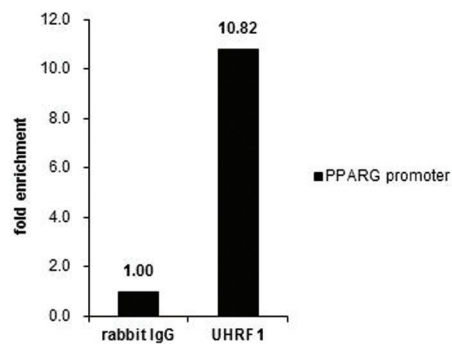
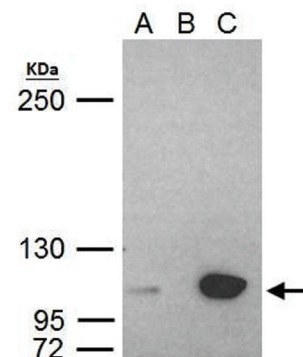


Figure 1. UHRF1 ChIP results

ChIP was performed with HCT116 chromatin extract and 5 µg of either control rabbit IgG or UHRF1 antibody. The precipitated DNA was detected by PCR with primer set targeting to PPARG promoter.

Figure 2. IP results

UHRF1 antibody immunoprecipitates UHRF1 protein in IP experiments. IP Sample: HeLa whole cell extract A. 40 µg HeLa whole cell extract B. Control with 2 µg of preimmune rabbit IgG C. Immunoprecipitation of UHRF1 protein by 2 µg of UHRF1 antibody (C15410258) 5% SDS-PAGE The immunoprecipitated UHRF1 protein was detected by western blot with the UHRF1 antibody (C15410258) diluted 1:1,000.



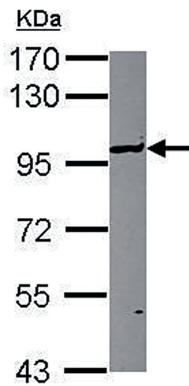


Figure 3. Western blot

Sample: 30 µg of HCT116 whole cell lysate
7.5% SDS PAGE
UHRF1 diluted 1:1,000

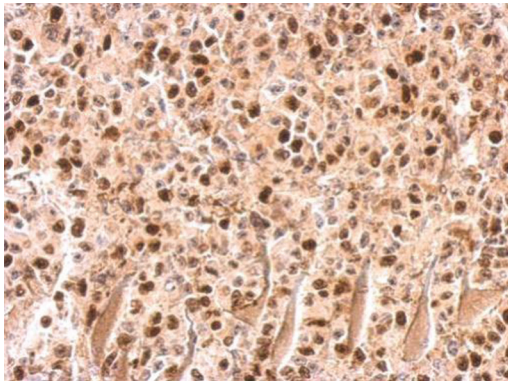


Figure 4. IHC

UHRF1 antibody detects UHRF1 protein on HBL435 xenograft by immunohistochemical analysis.
Sample: Paraffin-embedded HBL435 xenograft.
UHRF1 antibody (C15410258) dilution: 1:500.

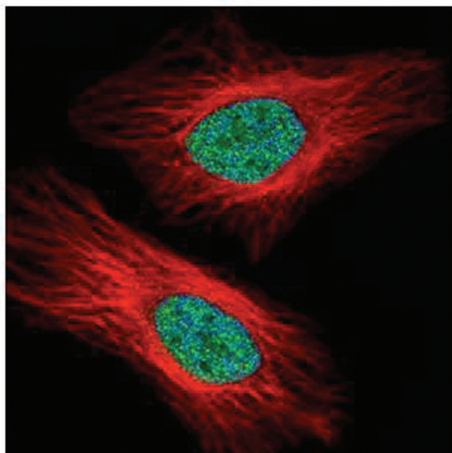


Figure 5. IFA

Confocal immunofluorescence analysis of paraformaldehyde-fixed HeLa cells using UHRF1 antibody (Cat. No. C15410258) (Green) at a 1:500 dilution. Red: Alpha-tubulin filaments.

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 20, 2021