

H4K8ac polyclonal antibody - Classic

Cat. No. C15410235

Type: Polyclonal

Source: Rabbit

Lot #: 41423

Size: 25 µl/100 µl

Concentration: 1 µg/µl

Specificity: Human, mouse, rat

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
Dot blot	1:5,000	Figure 3
ICC/IF	1:100-1:1,000	Figure 5
IHC	1:100-1:1,000	Figure 6, 7, 8
Immunoprecipitation	1:100-1:5,000	Figure 4
Western blot	1:1,000-1:10,000	Figure 2

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

Results

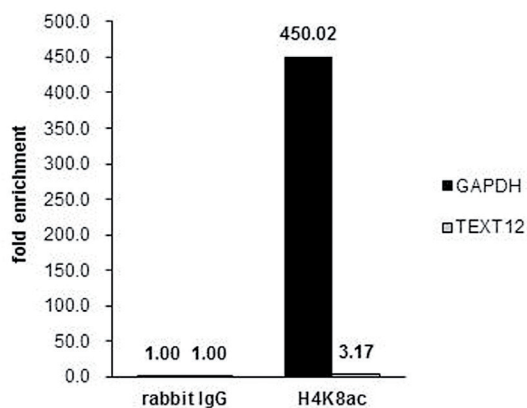


Figure 1. H4K8ac ChIP results

ChIP was performed with HeLa chromatin extract treated with Trichostatin A (0.4 µM for 18 h) and 5 µg of either control rabbit IgG or H4K8ac antibody. The precipitated DNA was detected by PCR with primer set targeting to GAPDH or TEXT12.

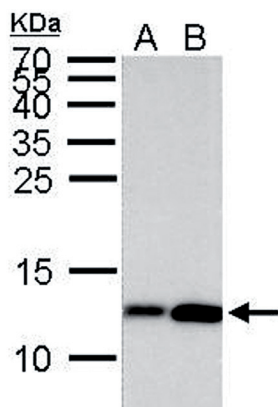


Figure 2. H4K8ac WB results

H4K8ac antibody detects Histone H4 acetylated at Lys8 by Western blot analysis.

A. 30 μg HeLa whole extract (0.012% DMSO treatment for 18 hr)

B. 30 μg HeLa whole extract (0.4 μM Trichostatin A treatment for 18 hr)

15 % SDS-PAGE

H4K8ac antibody dilution: 1:5,000

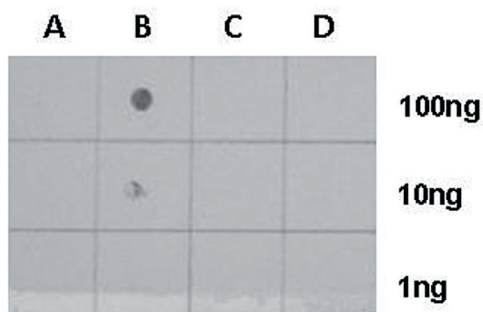


Figure 3. H4K8ac Dot blot results

Dotblot analysis of H4K8ac antibody with different peptides.

Varied amount of the peptides were spotted onto positively charged nylon membrane and blotted with the H4K8ac antibody at a 1:5,000 dilution.

A: Peptide H4K8un

B: Peptide H4K8ac

C: Peptide H4K16un

D: Peptide H4K16ac

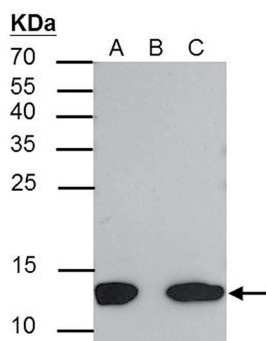


Figure 4. H4K8ac IP results

H4K8ac antibody immunoprecipitates Histone H4 acetylated at Lys8 in IP experiments.

IP Sample: HeLa whole cell extract A: 30 μg whole cell extract of HeLa cells (0.4 μM TSA treatment for 18hr) B: Control with 2.5 μg of pre-immune rabbit IgG C: Immunoprecipitation of Histone H4 acetylated at Lys8 by 2.5 μg of H4K8ac antibody. The immunoprecipitated Histone H4 (acetyl K8) protein was detected by western blot with the H4K8ac antibody diluted 1:5,000.

TECHNICAL DATASHEET

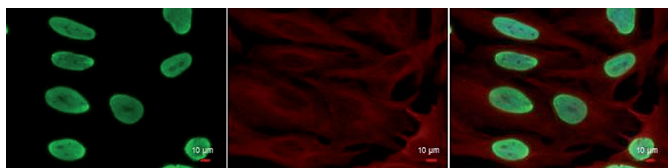


Figure 5. H4K8ac IF results

H4K8ac antibody detects Histone H4 acetylated at Lys8 in the nucleus by immunofluorescent analysis.

Sample: HeLa cells were fixed in 4% paraformaldehyde at RT for 15 min.

Green: Histone H4 acetylated at Lys8 stained by H4K8ac antibody diluted 1:500.

Red: alpha Tubulin antibody, a cytoskeleton marker.

Blue: Hoechst 33342 staining.

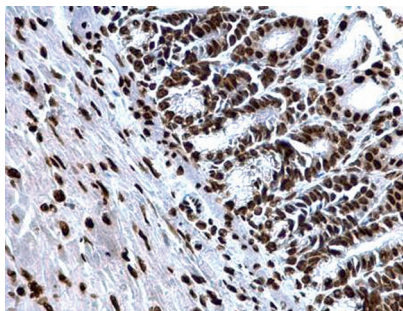


Figure 6. H4K8ac IHC results

H4K8ac antibody detects Histone H4 acetylated at Lys8 at nucleus on mouse duodenum by immunohistochemical analysis.

Sample: Paraffin-embedded mouse duodenum.

H4K8ac antibody dilution: 1:500.

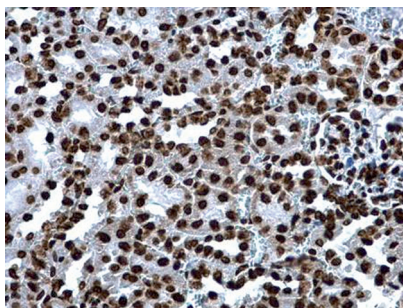


Figure 7. H4K8ac IHC results

H4K8ac antibody detects Histone H4 acetylated at Lys8 at nucleus on mouse kidney by immunohistochemical analysis.

Sample: Paraffin-embedded mouse kidney.

H4K8ac antibody dilution: 1:500.

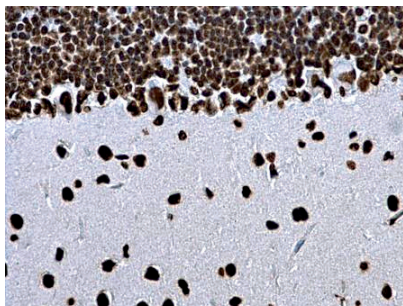


Figure 8. H4K8ac IHC results

H4K8ac antibody detects Histone H4 acetylated at Lys8 at nucleus on rat hind brain by immunohistochemical analysis.

Sample: Paraffin-embedded rat hind brain.

H4K8ac antibody dilution: 1:500.

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