

## H3K27ac antibody

**Cat. No.** C15210016

**Type:** Monoclonal, ChIP grade/ChIP-seq grade

**Source:** Rabbit

**Lot:** 006

**Size:** 100 µg

**Concentration:** 1 µg/µl

**Specificity:** Human, wide range expected

**Purity:** Affinity purified antibody.

**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

**Storage buffer:** PBS containing 50% glycerol, 1% BSA and 0.09% azide

**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Description:** Monoclonal antibody raised in rabbit against the region of histone H3 containing the acetylated lysine 27 (H3K27ac), using a KLH-conjugated synthetic peptide.

### Applications

Applications	Suggested dilution	References
ChIP/ChIP-seq*	0.5 - 1 µg per IP	Fig 1, 2
Dot Blotting	1:2,000	Fig 3
Western Blotting	1:1,000	Fig 4
Immunofluorescence	1:1,000	Fig 5

\*Please note that the optimal antibody amount per IP should be determined by the end-user. We recommend testing 0.5 - 5 µg per IP.

### Target description

Histones are the main constituents of the protein part of chromosomes of eukaryotic cells. They are rich in the amino acids arginine and lysine and have been greatly conserved during evolution. Histones pack the DNA into tight masses of chromatin. Two core histones of each class H2A, H2B, H3 and H4 assemble and are wrapped by 146 base pairs of DNA to form one octameric nucleosome. Histone tails undergo numerous post-translational modifications, which either directly or indirectly alter chromatin structure to facilitate transcriptional activation or repression or other nuclear processes. In addition to the genetic code, combinations of the different histone modifications reveal the so-called "histone code". Histone methylation and demethylation is dynamically regulated by respectively histone methyl transferases and histone demethylases. Acetylation of histone H3K27 is associated with active genes.

#### Diagenode sa. BELGIUM | EUROPE

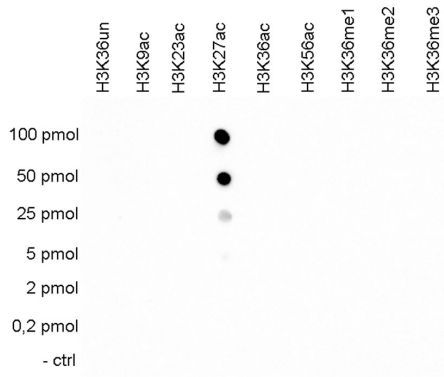
LIEGE SCIENCE PARK  
Rue du Bois Saint-Jean, 3  
4102 Seraing - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

#### Diagenode LLC. 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

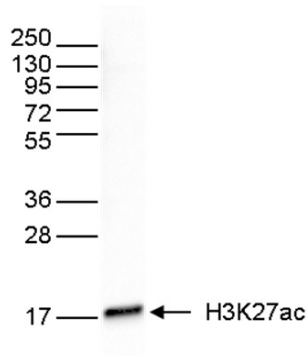
Last update: January, 2024





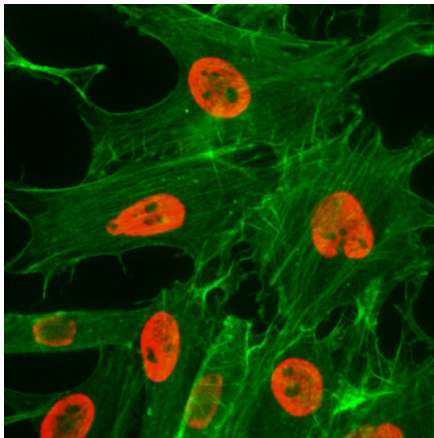
**Figure 3. Cross reactivity tests using the Diagenode monoclonal antibody directed against H3K27ac**

To test the cross reactivity of the Diagenode antibody against H3K27ac (cat. No. C15210016), a Dot Blot analysis was performed with peptides containing other histone modifications and the unmodified H3K27. One hundred to 0.2 pmol of the respective peptides were spotted on a membrane. The antibody was used at a dilution of 1:2,000. Figure 3 shows a high specificity of the antibody for the modification of interest.



**Figure 4. Western blot analysis using the Diagenode monoclonal antibody directed against H3K27ac**

Western blot was performed on whole cell extracts (40 µg) from HeLa cells using the Diagenode antibody against H3K27ac (cat. No. C15210016). The antibody was diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is shown on the right, the marker (in kDa) is shown on the left.



**Figure 5. Immunofluorescence using the Diagenode monoclonal antibody directed against H3K27ac**

HeLa cells were stained with the Diagenode antibody against H3K27ac (cat. No. C15210016, red) diluted 1:1,000. Actin was stained with fluorescein phalloidin (green).