

## H3K56ac antibody

**Cat. No.** **C15410213**

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

**Source:** Rabbit

**Lot:** A2069P

**Size:** 50 µg

**Concentration:** 0.8 µg/µl

**Specificity:** Human: positive  
Other species: not tested.

**Purity:** Affinity purified polyclonal antibody.

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

**Storage buffer:** PBS containing 0.05% azide and 0.05% ProClin 300.

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

**Description:** Polyclonal antibody raised in rabbit against the region of histone H3 containing the acetylated lysine 56 (H3K56ac), using a KLH-conjugated synthetic peptide.

### Applications

Applications	Suggested dilution	References
ChIP*	1 µg per ChIP	Fig 1, 2
CUT&TAG	1 µg	Fig 3
ELISA	1:1000	Fig 4
Dot Blotting	1:20,000	Fig 5
Western Blotting	1:200	Fig 6

\*Please note that the optimal antibody amount per IP should be determined by the end-user. We recommend testing 1-5 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.

#### Diagenode sa. BELGIUM | EUROPE

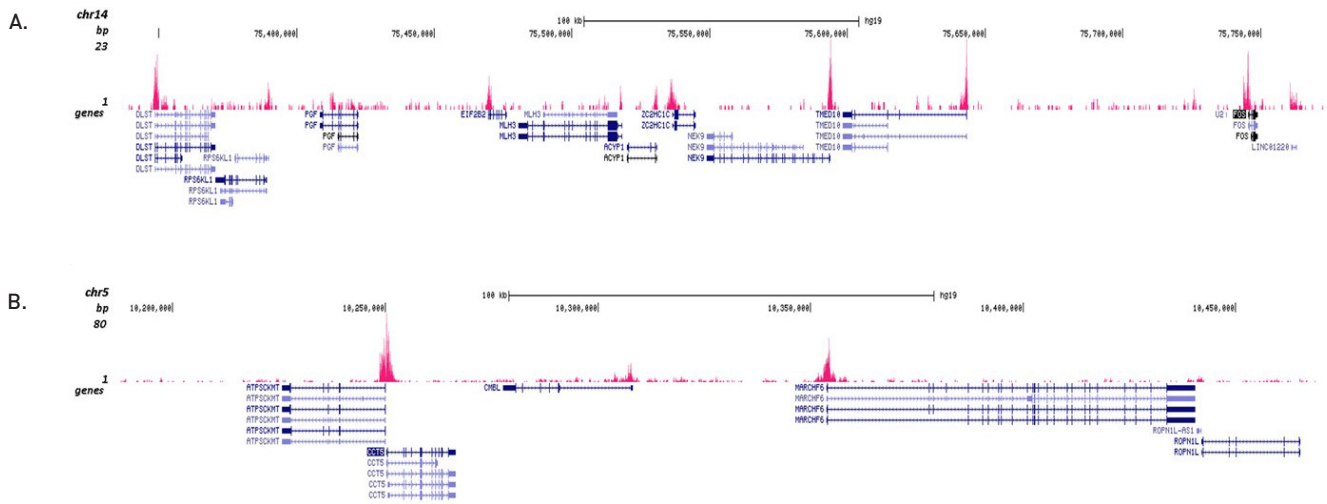
LIEGE SCIENCE PARK  
Rue du Bois Saint-Jean, 3  
4102 Seraing - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

#### 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

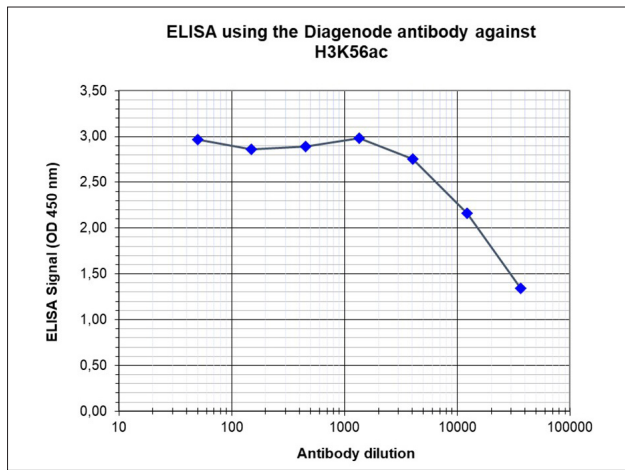
Last update: February, 2023





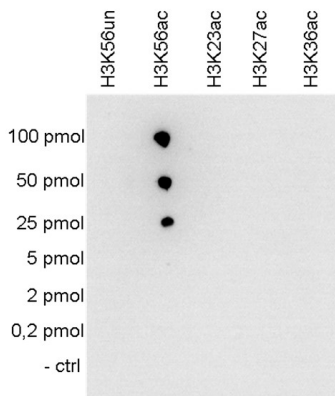
**Figure 3. Cut&Tag results obtained with the Diagenode antibody directed against H3K56ac**

CUT&TAG (Kaya-Okur, H.S., Nat Commun 10, 1930, 2019) was performed on 50,000 K562 cells using 1 µg of the Diagenode antibody against H3K56ac (cat. No. C15410213) and the Diagenode pA-Tn5 transposase (C01070001). The libraries were subsequently analysed on an Illumina NextSeq 500 sequencer (2x75 paired-end reads) according to the manufacturer’s instructions. The tags were aligned to the human genome (hg19) using the BWA algorithm. Figure 3 shows the peak distribution in 2 genomic regions surrounding the FOS gene on chromosome 14 and the CCT5 gene on chromosome 5 (figure 3A and B, respectively).



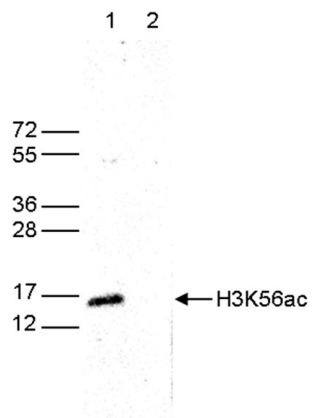
**Figure 4. Determination of the antibody titer**

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against H3K56ac (cat. No. 15410213) in antigen coated wells. The antigen used was a peptide containing the histone modification of interest. By plotting the absorbance against the antibody dilution (Figure 4), the titer of the antibody was estimated to be 1:36,500.



**Figure 5. Cross reactivity tests using the Diagenode antibody directed against H3K56ac**

To test the cross reactivity of the Diagenode antibody against H3K56ac (cat. No. 15410213), a Dot Blot analysis was performed with peptides containing other histone modifications and the unmodified H3K56. One hundred to 0.2 pmol of the respective peptides were spotted on a membrane. The antibody was used at a dilution of 1:20,000. Figure 5 shows a high specificity of the antibody for the modification of interest.



**Figure 6. Western blot analysis using the Diagenode antibody directed against H3K56ac**

Western blot was performed on whole cell extracts (40 µg, lane 1) from HeLa cells, and on 1 µg of recombinant histone H3 (lane 2) using the Diagenode antibody against H3K56ac (cat. No. C15410213). The antibody was diluted 1:200 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.