

## H3K79ac monoclonal antibody

**Cat. No.** C15210006

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

**Source:** Rabbit

**Lot #:** 001

**Size:** 100 µg/100 µl

**Concentration:** 1 µg/µl

**Specificity:** Human: positive

Other species: not tested

**Purity:** Protein A purified monoclonal antibody in PBS containing 1% BSA and 0.09% azide.

**Storage:** Store at -4°C.

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

**Description:** Monoclonal antibody raised in rabbit against histone H3 acetylated at Lys79 (H3K79ac), using a KLH-conjugated synthetic peptide.

### Applications

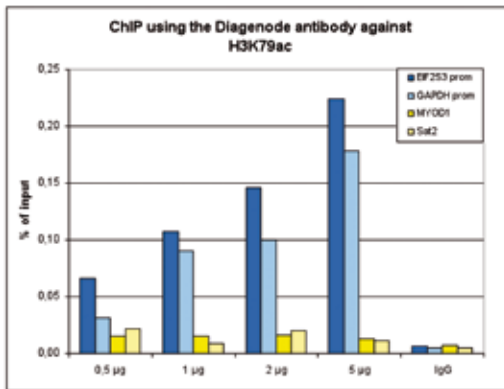
	Suggested dilution	Results
ChIP*	5 µg/ChIP	Fig 1, 2
Western blotting	1:1,000	Fig 3
Immunofluorescence	1:500	Fig 4

\* Please note that the optimal antibody amount per IP should be determined by the end-user. We recommend testing 1-10 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 Lys79 is associated with gene activation.

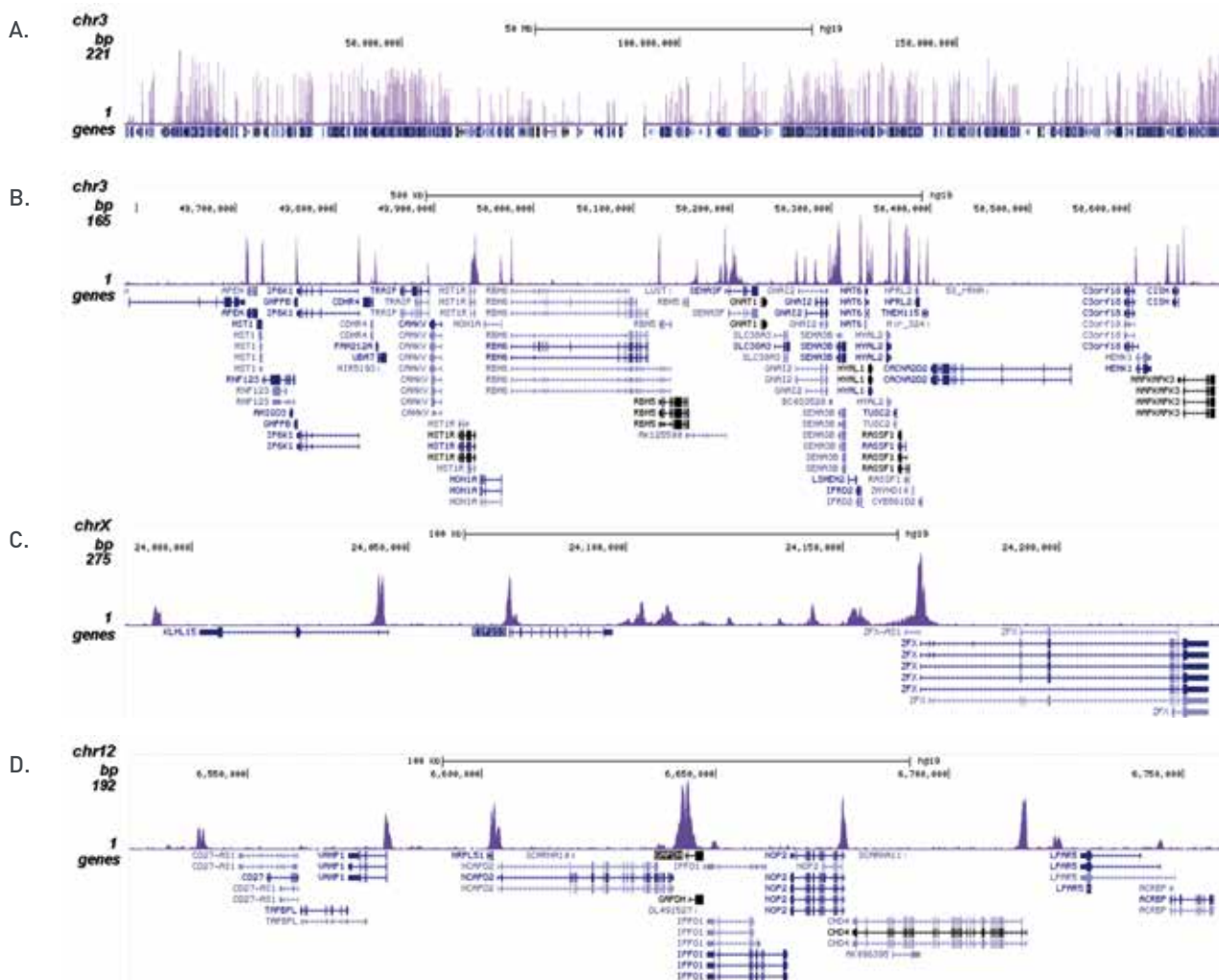
## Results



**Figure 1. ChIP results obtained with the Diagenode monoclonal antibody directed against H3K79ac**

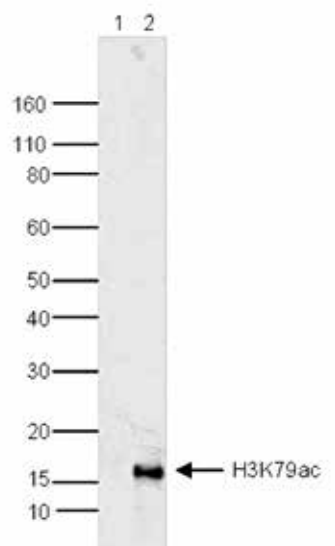
ChIP assays were performed using HeLa cells, the Diagenode antibody against H3K79ac [cat. No. C15210006] and optimized PCR primer sets for qPCR. ChIP was performed with the “iDeal ChIP-seq” kit [cat. No. C01010051], using sheared chromatin from 1 million cells. A titration consisting of 0.5, 1, 2 and 5 µg of antibody per ChIP experiment was analyzed. IgG (1 µg/IP) was used as a negative IP control. Quantitative PCR was performed with optimized primers for the promoters of the EIF2S3 and GAPDH genes, used as positive control, and for the MYOD1 gene and the Sat2 satellite repeat, used as negative control.

Figure 1 shows the recovery, expressed as a % of input (the relative amount of immunoprecipitated DNA compared to input DNA after qPCR analysis).



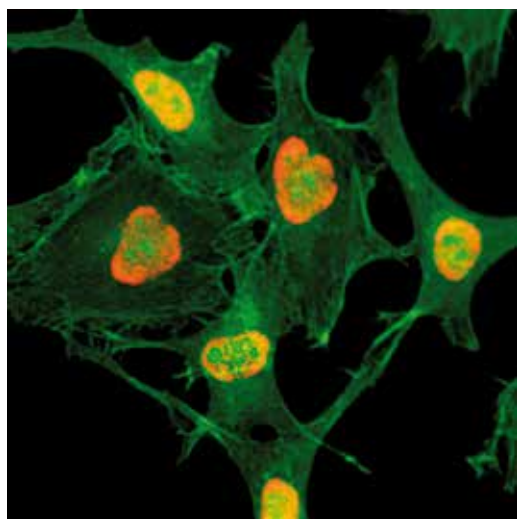
**Figure 2. ChIP-seq results obtained with the Diagenode monoclonal antibody directed against H3K79ac**

ChIP was performed on sheared chromatin from 1 million HeLa cells using 5 µg of the Diagenode antibody against H3K79ac (Cat. No. C15210006) as described above. The IP'd DNA was subsequently analysed on an Illumina HiSeq. Library preparation, cluster generation and sequencing were performed according to the manufacturer's instructions. The 50 bp tags were aligned to the human genome using the BWA algorithm. Figure 2 shows the enrichment along the complete sequence and a 1 Mb region of human chromosome 3 (fig 2A and B), and in two genomic regions surrounding the EIF2S3 and GAPDH positive control genes (fig 2C and D).



**Figure 3. Western blot analysis using the Diagenode monoclonal antibody directed against H3K79ac**

Histone extracts from HeLa cells treated with butyrate (lane 2) or untreated control cells (lane 1) were analysed by Western blot using the Diagenode monoclonal antibody against H3K79ac (Cat. No. C15210006) diluted 1:1,000 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.



**Figure 4. Immunofluorescence using the Diagenode monoclonal antibody directed against H3K79ac**

HeLa cells treated with butyrate were stained with the Diagenode antibody against H3K79ac (Cat. No. C15210006 red) diluted 1:500. Actin filaments were stained with fluorescein phalloidin (green).

**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 20, 2016