



H3K4me1 polyclonal antibody

Cat. No. C15410194 (pAb-194-050)

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

Source: Rabbit

Lot #: A1862D

Size: 50 µg/ 34 µl

Concentration: 1.5 µg/µl

Specificity: Human, mouse, wide range: expected

Purity: Affinity purified polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

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.

Description: Polyclonal antibody raised in rabbit against the region of histone H3 containing the monomethylated lysine 4 (H3K4me1), using a KLH-conjugated synthetic peptide.

Applications

	Suggested dilution	Results
ChIP*	0.5 - 1 µg per IP	Fig 1, 2
ELISA	1:400	Fig 3
Dot blotting/Peptide array	1:5,000/1:2,000	Fig 4
Western blotting	1:500	Fig 5
IF	1:200	Fig 6

* 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. Methylation of histone H3K4 is associated with active genes.

Results

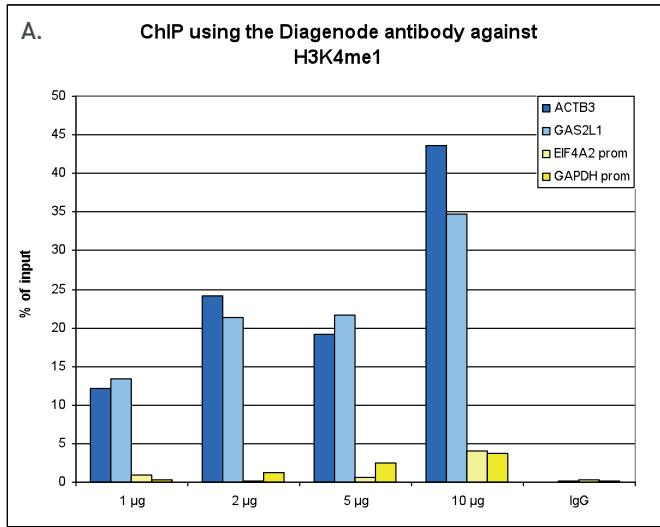


Figure 1. ChIP results obtained with the Diagenode antibody directed against H3K4me1

ChIP assays were performed using human K562 cells, the Diagenode antibody against H3K4me1 (Cat. No. C15410194) and optimized PCR primer pairs for qPCR. ChIP was performed with the “iDeal ChIP-seq” kit (Cat. No. AB-001-0024), using sheared chromatin from 100,000 cells. A titration consisting of 1, 2, 5 and 10 µg of antibody per ChIP experiment was analyzed. IgG (1 µg/IP) was used as a negative IP control. Quantitative PCR was performed with primers for a region surrounding the ACTB and GAS2L1 gene, respectively, used as positive controls, and for the promoters of the GAPDH and EIF4A2 genes, used as negative controls. Figure 1 shows the recovery, expressed as a % of input (the relative amount of immunoprecipitated DNA compared to input DNA after qPCR analysis).

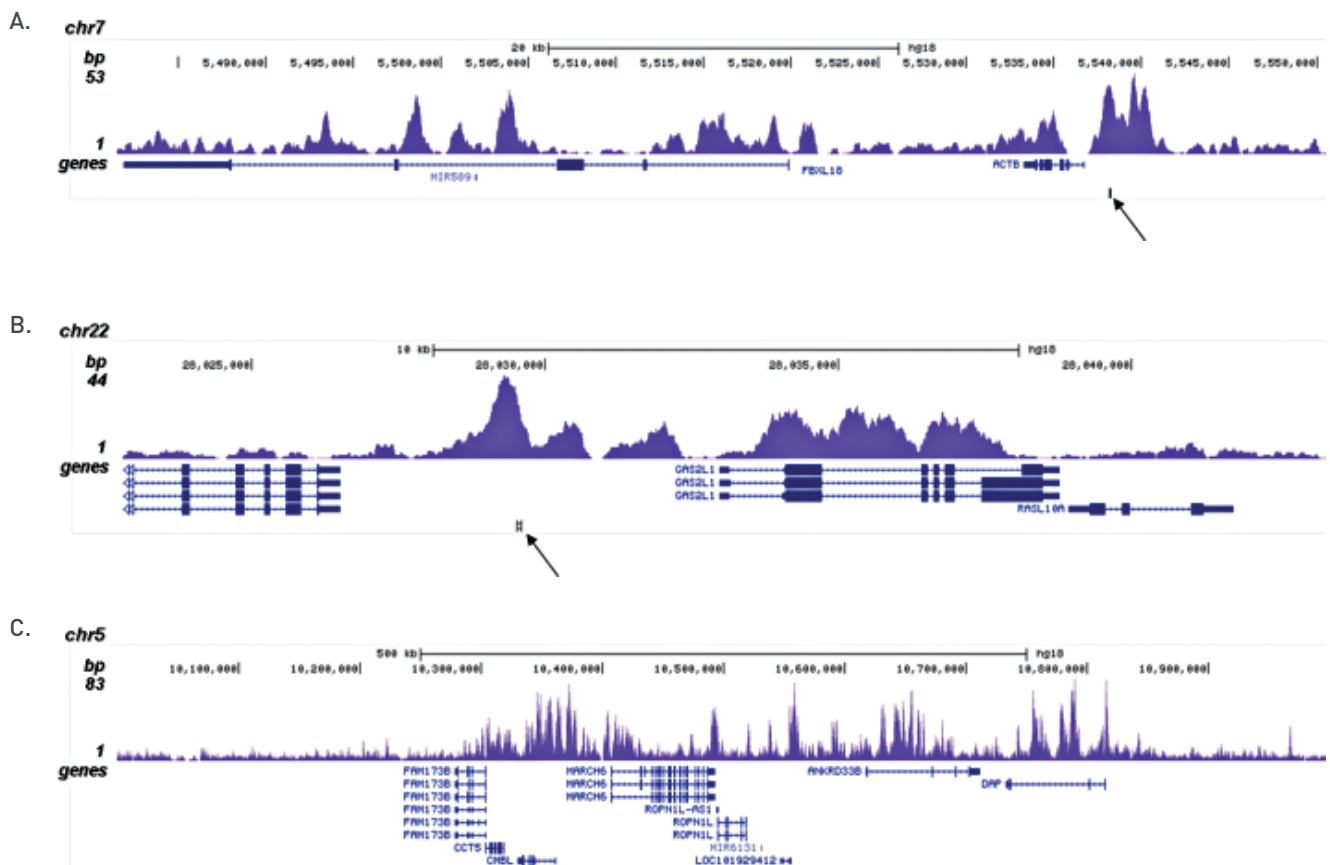


Figure 2. ChIP-seq results obtained with the Diagenode antibody directed against H3K4me1

ChIP was performed on sheared chromatin from 100,000 K562 cells with the “iDeal ChIP-seq” kit (Cat. No. AB-001-0024) using 1 µg of the Diagenode antibody against H3K4me1 (Cat. No. C15410194) as described above. The IP’d DNA was subsequently analysed on an Illumina Genome Analyzer. Library preparation, cluster generation and sequencing were performed according to the manufacturer’s instructions. The 36 bp tags were aligned to the human genome using the ELAND algorithm. Figure 2A and B show the H3K4me1 signal in two genomic regions containing the ACTB and GAS2L1 positive controls. The position of the amplicon used for ChIP-qPCR is indicated by an arrow. Figure 2C shows the H3K4me1 peak distribution along a 1 Mb genomic region of chromosome 5.

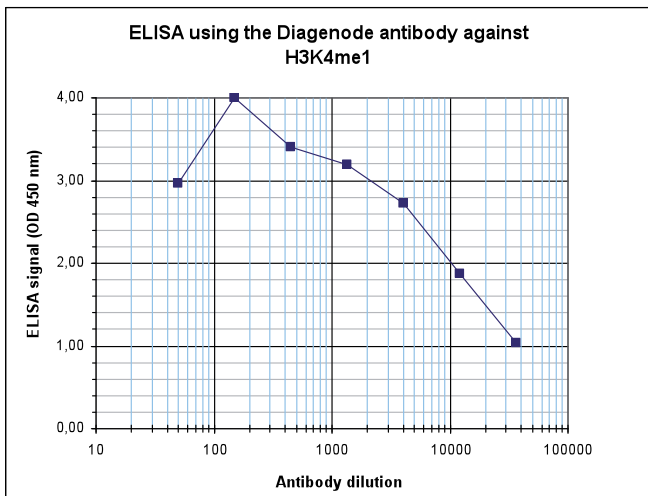


Figure 3. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody against H3K4me1 [Cat. No. C15410194]. The antigen used was a peptide containing the histone modification of interest. By plotting the absorbance against the antibody dilution (Figure 3), the titer of the antibody was estimated to be 1:10,300.

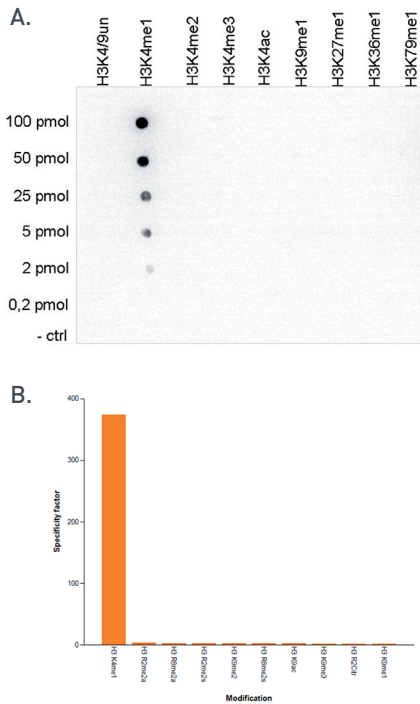


Figure 4. Cross reactivity tests using the Diagenode antibody directed against H3K4me1

Figure 4A To test the cross reactivity of the Diagenode antibody against H3K4me1 [Cat. No. C15410194], a Dot Blot analysis was performed with peptides containing other histone modifications and the unmodified H3K4. One hundred to 0.2 pmol of the respective peptides were spotted on a membrane. The antibody was used at a dilution of 1:5,000. Figure 4A shows a high specificity of the antibody for the modification of interest.

Figure 4B The specificity of the antibody was further demonstrated by peptide array analyses on an array containing 384 peptides with different combinations of modifications from histone H3, H4, H2A and H2B. The antibody was used at a dilution of 1:2,000. Figure 4B shows the specificity factor, calculated as the ratio of the average intensity of all spots containing the mark, divided by the average intensity of all spots not containing the mark.

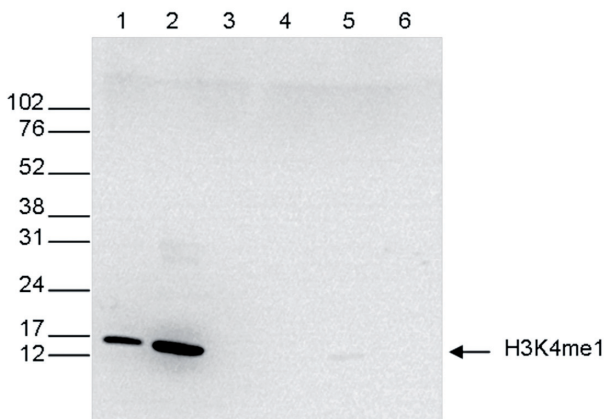


Figure 5. Western blot analysis using the Diagenode antibody directed against H3K4me1

Western blot was performed on whole cell (25 µg, lane 1) and histone extracts (15 µg, lane 2) from HeLa cells, and on 1 µg of recombinant histone H2A, H2B, H3 and H4 (lane 3, 4, 5 and 6, respectively) using the Diagenode antibody against H3K4me1 [Cat. No. C15410194]. The antibody was diluted 1:500 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.

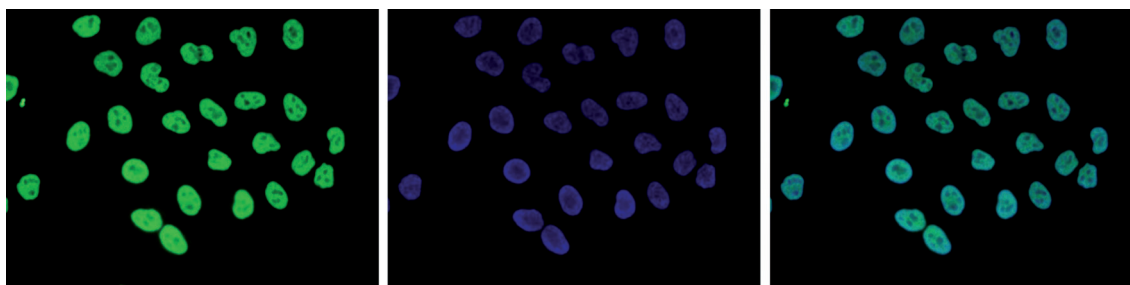


Figure 6. Immunofluorescence using the Diagenode antibody directed against H3K4me1

HeLa cells were stained with the Diagenode antibody against H3K4me1 (Cat. No. C15410194) and with DAPI. Cells were fixed with 4% formaldehyde for 10' and blocked with PBS/TX-100 containing 5% normal goat serum and 1% BSA. The cells were immunofluorescently labeled with the H3K4me1 antibody (left) diluted 1:200 in blocking solution followed by an anti-rabbit antibody conjugated to Alexa488. The middle panel shows staining of the nuclei with DAPI. A merge of the two stainings is shown on the right.

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: March 18, 2014