

## Histone H3K18ac (Acetyl H3K18) Polyclonal Antibody

(Catalog # A-4024)

### Background

Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. Nucleosomes consist of approximately 146 bp of DNA wrapped around a histone octamer composed of pairs of each of the four core histones (H2A, H2B, H3, and H4). The chromatin fiber is further compacted through the interaction of a linker histone, H1, with the DNA between the nucleosomes to form higher order chromatin structures. This gene is intronless and encodes a member of the histone H3 family. Transcripts from this gene lack polyA tails; instead, they contain a palindromic termination element. This gene is located separately from the other H3 genes that are in the histone gene cluster on chromosome 6p22-p21. 3.

### Description

Histone H3K18ac (Acetyl H3K18) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.05% proclin300, 50% glycerol, pH7.3

### Specificity

Broad Range, Mouse, Rat, Human

### Isotype

IgG

### Purification

Affinity Purified

### Immunogen

A synthetic acetylated peptide around K9 of human Histone H3 (NP\_003520.1).

### Storage

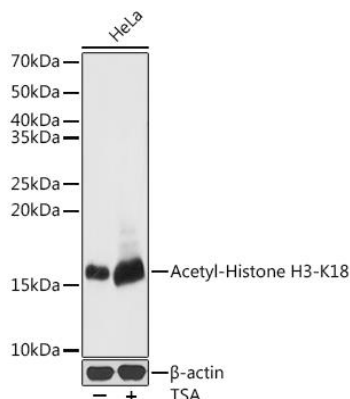
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

### Alternative Names

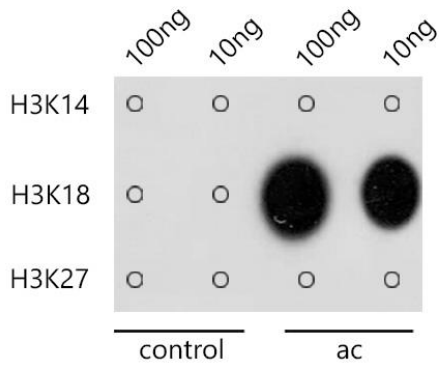
H3K18ac antibody; H3K18a antibody; H3t; H3.4; H3/g; H3FT

### Application

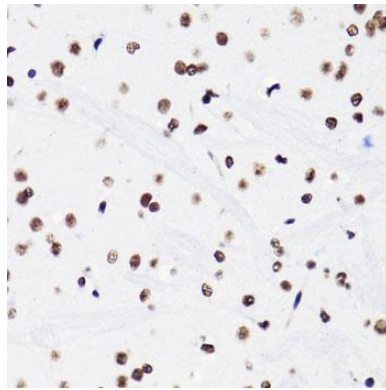
WB, IHC, IF/ICC, IP, ChIP, ChIPseq; Recommended dilution: WB 1:100 - 1:500, IHC 1:200 - 1:500, IF/ICC 1:500 - 1:1000, IP 1:200 - 1:500, ChIP 1:50 - 1:100, ChIP-seq 1:50 - 1:100



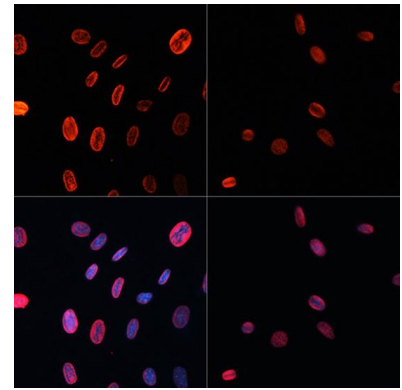
Western blot analysis of extracts of HeLa cells, using Acetyl-Histone H3-K18 antibody at 1:500 dilution. HeLa cells were treated by TSA (1  $\mu$ M) at 37°C for 18 hours. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25 $\mu$ g per lane. Blocking buffer: 3% nonfat dry milk in TBST.



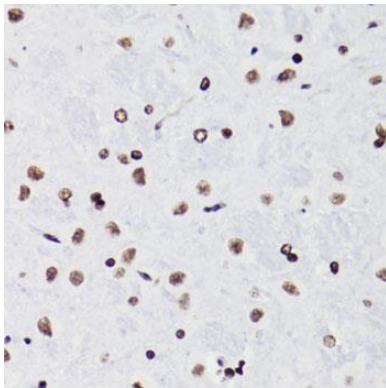
Dot-blot analysis of all sorts of methylation peptides using Acetyl-Histone H3-K18 antibody at 1:1000 dilution.



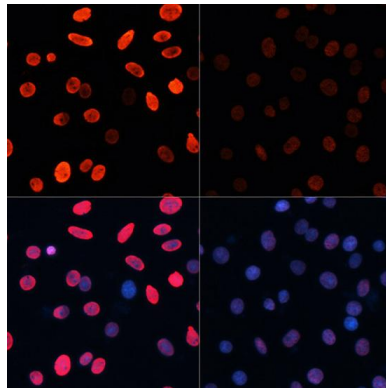
Immunohistochemistry of paraffin-embedded mouse brain using H3K18ac antibody at dilution of 1:100 (40x lens).



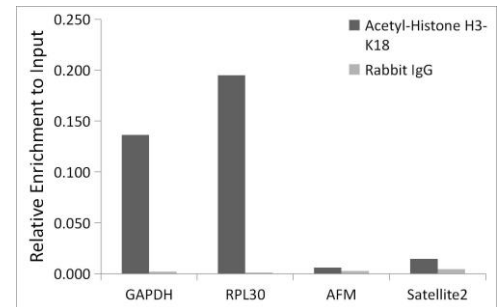
Immunofluorescence analysis of NIH/3T3 cells using Acetyl-Histone H3-K18 antibody at dilution of 1:100. NIH/3T3 cells were treated by TSA (1  $\mu$ M) at 37 degrees Celsius for 18 hours. Blue: DAPI for nuclear staining.



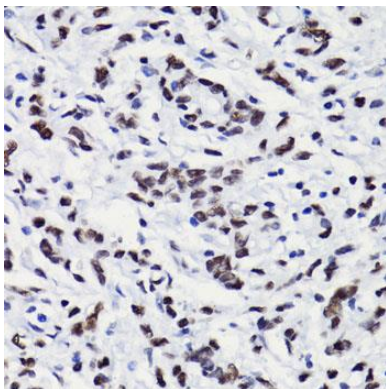
Immunohistochemistry of paraffin-embedded rat brain using H3K18ac antibody at dilution of 1:100 (40x lens).



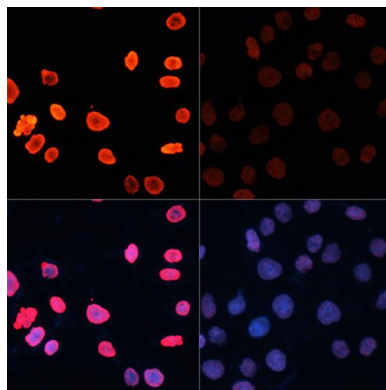
Immunofluorescence analysis of C6 cells using Acetyl-Histone H3-K18 antibody at dilution of 1:100. C6 cells were treated by TSA (1  $\mu$ M) at 37 degrees Celsius for 18 hours. Blue: DAPI for nuclear staining.



Chromatin immunoprecipitation analysis of extracts of HCT116 cells, using Acetyl-Histone H3-K18 Rabbit pAb antibody and rabbit IgG. The amount of immunoprecipitated DNA was checked by quantitative PCR. Histogram was constructed by the ratios of the immunoprecipitated DNA to the input.



Immunohistochemistry of paraffin-embedded human gastric cancer using H3K18ac antibody at dilution of 1:100 (40x lens).



Immunofluorescence analysis of HeLa cells using Acetyl-Histone H3-K18 antibody at dilution of 1:100. HeLa cells were treated by TSA (1  $\mu$ M) at 37 degrees Celsius for 18 hours. Blue: DAPI for nuclear staining.