
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

Uniprot ID

Q16695/P68431

Purification

Affinity Purified

Immunogen

A synthetic acetylated peptide around K18 of human H3 (NP_003520.1).

Storage

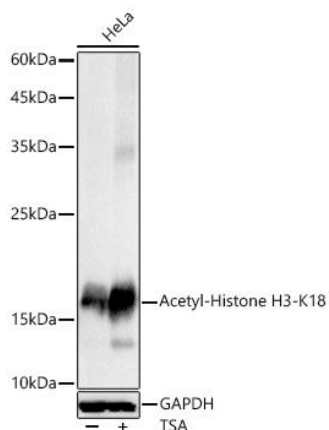
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

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

Application

WB, IHC, IF/ICC, ChIP, DB, ELISA; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:200, IF/ICC 1:50 - 1:200, ChIP 5µg antibody for 5µg-10µg of Chromatin, DB 1:500 - 1:1000, ELISA - recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.



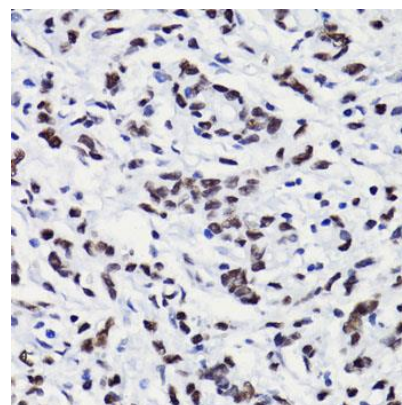
Western blot analysis of lysates from HeLa cells using Histone H3K18ac (Acetyl H3K18) Polyclonal Antibody at 1:1000 dilution. HeLa cells were treated by TSA (1 μ M) at 37°C for 18 hours.

Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.

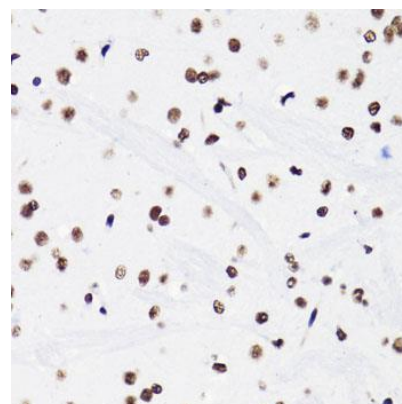
Lysates/proteins: 25 μ g per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

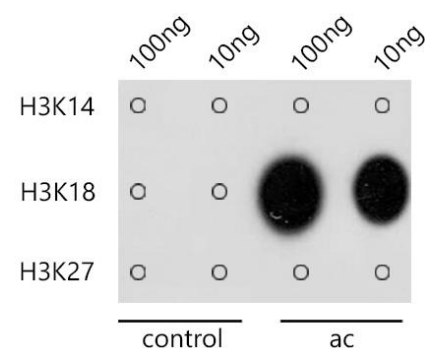
Exposure time: 0.5s.



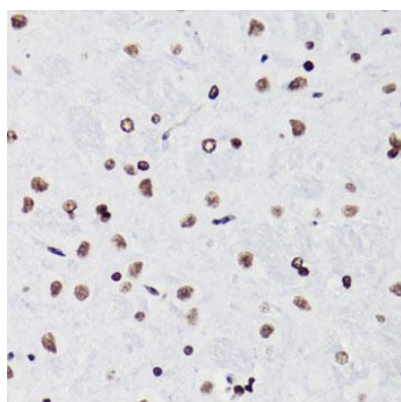
Immunohistochemistry analysis of paraffin-embedded Human gastric cancer using Histone H3K18ac (Acetyl H3K18) Polyclonal Antibody at dilution of 1:100 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



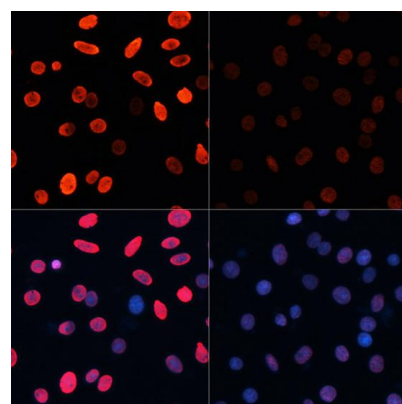
Immunohistochemistry analysis of paraffin-embedded Mouse brain using Histone H3K18ac (Acetyl H3K18) Polyclonal Antibody at dilution of 1:100 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



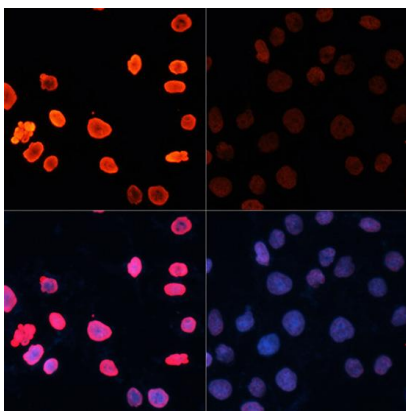
Dot-blot analysis of all sorts of methylation peptides using Histone H3K18ac (Acetyl H3K18) Polyclonal Antibody at 1:1000 dilution.



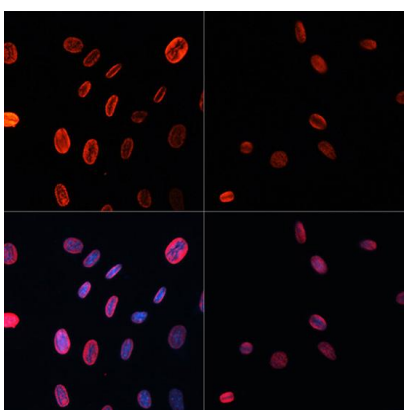
Immunohistochemistry analysis of paraffin-embedded Rat brain using Histone H3K18ac (Acetyl H3K18) Polyclonal Antibody at dilution of 1:100 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



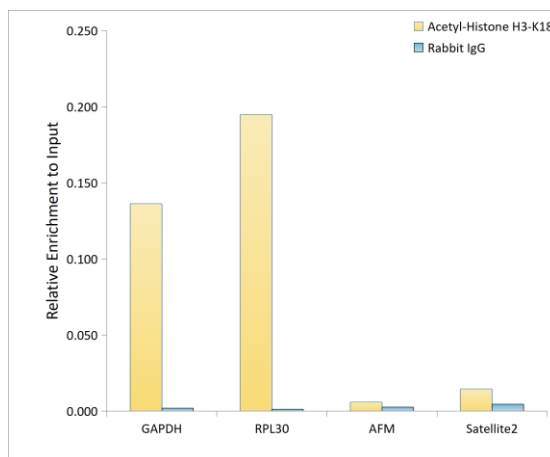
Immunofluorescence analysis of C6 cells using Histone H3K18ac (Acetyl H3K18) Polyclonal Antibody at dilution of 1:100. C6 cells were treated by TSA (1 μ M) at 37°C for 18 hours. Blue: DAPI for nuclear staining.



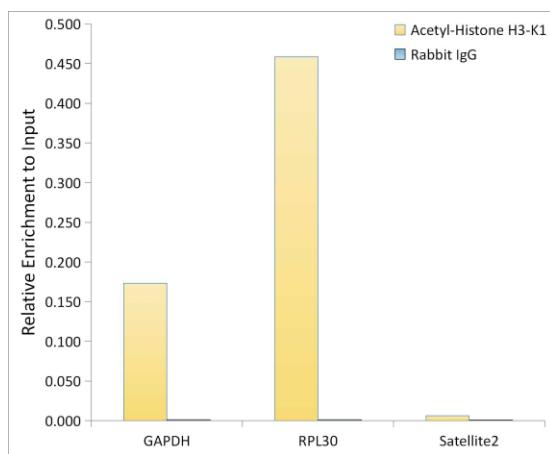
Immunofluorescence analysis of HeLa cells using Histone H3K18ac (Acetyl H3K18) Polyclonal Antibody at dilution of 1:100. HeLa cells were treated by TSA (1 μ M) at 37°C for 18 hours. Blue: DAPI for nuclear staining.



Immunofluorescence analysis of NIH/3T3 cells using Histone H3K18ac (Acetyl H3K18) Polyclonal Antibody at dilution of 1:100. NIH/3T3 cells were treated by TSA (1 μ M) at 37°C for 18 hours. Blue: DAPI for nuclear staining.



Chromatin immunoprecipitation analysis of extracts of HCT116 cells, using Histone H3K18ac (Acetyl H3K18) Polyclonal 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.



Chromatin immunoprecipitation analysis of extracts of HCT116 cells, using Histone H3K18ac (Acetyl H3K18) Polyclonal 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.