
Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody

(Catalog # A-4022)

Background

Histone H3- along with H2A, H2B and H4- is involved in the structure of chromatin in eukaryotic cells. Histone H3 can undergo several different types of epigenetic modifications that influence cellular processes. These modifications including acetylation, phosphorylation, methylation, ubiquitination, and ADP-ribosylation occur on the N-terminal tail domains of histone H3, which results in remodeling of the nucleosome structure into an open conformation more accessible to transcription complexes. In most species, histone H3 is primarily acetylated at lysine 9, 14, 18, and 23.

Description

Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation:

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH 7.3.

Specificity

Broad Range, Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q16695/P68431

Purification

Affinity Purified

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 1-100 of human Histone H3 (NP_003520.1).

Storage

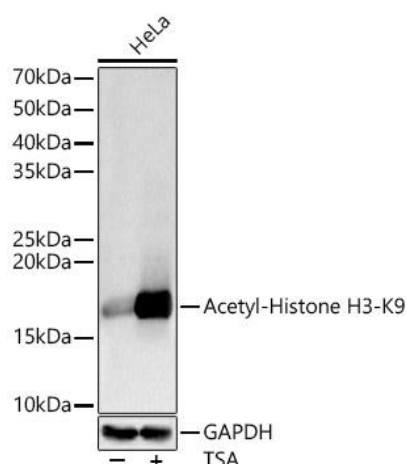
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

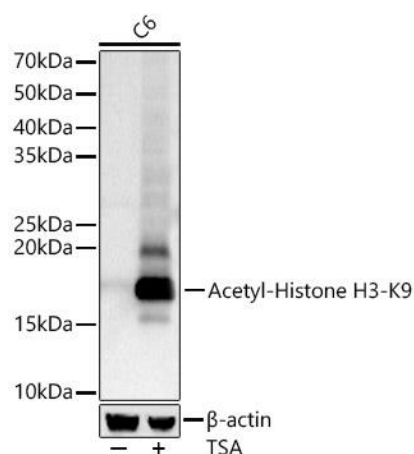
H3K9ac antibody, H3K9a antibody

Application

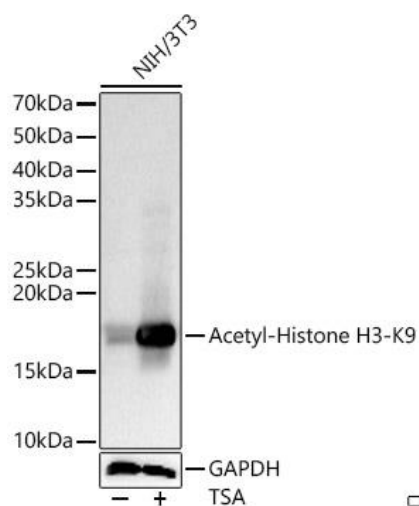
WB, IHC, IF/ICC, IP, ChIP, ChIP-seq, ELISA; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:200, IF/ICC 1:50 - 1:200, IP 0.5ug-4ug antibody for 200ug-400ug extracts of whole cells, ChIP 5µg antibody for 5µg-10µg of Chromatin, CHIP-seq 1:20 - 1:50, ELISA - Recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.



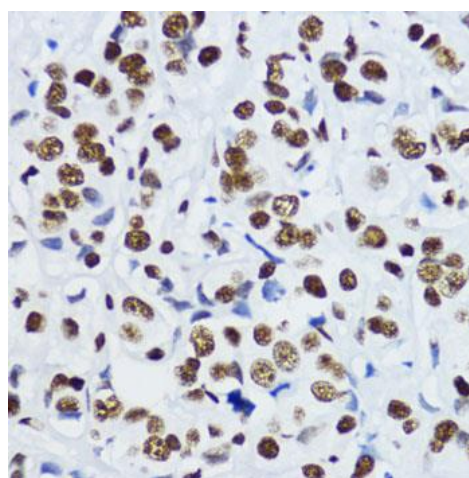
Western blot analysis of lysates from HeLa cells, using Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody at 1:1000 dilution. HeLa cells were treated by TSA (1 μ M) at 37°C for 18 hours. Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25 μ g per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 10s.



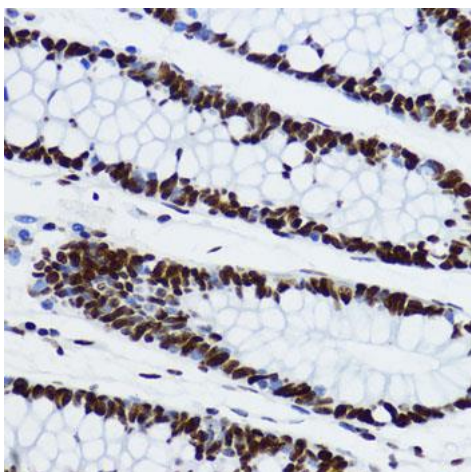
Western blot analysis of lysates from C6 cells, using Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody at 1:1000 dilution. C6 cells were treated by TSA (1 μ M) at 37°C for 18 hours. Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25 μ g per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 10s.



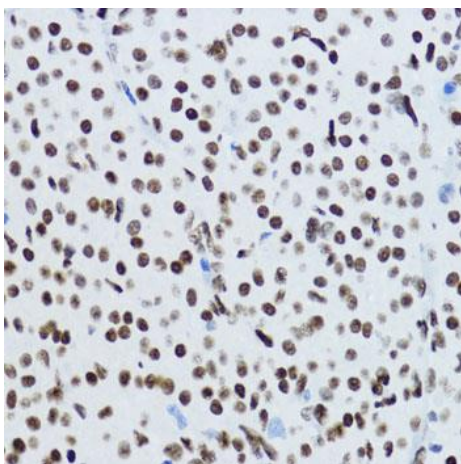
Western blot analysis of lysates from NIH/3T3 cells, using Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody at 1:1000 dilution. NIH/3T3 cells were treated by TSA (1 μ M) at 37°C for 18 hours. Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25 μ g per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 10s.



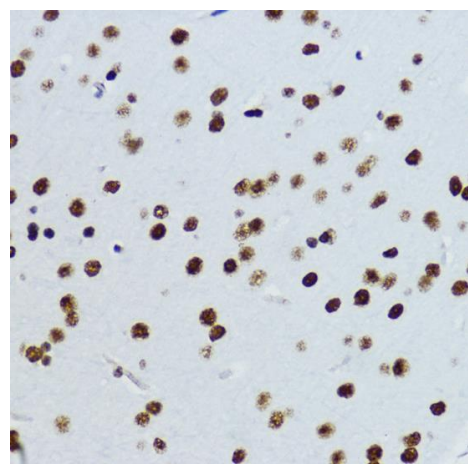
Immunohistochemistry analysis of paraffin-embedded Human mammary cancer using Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



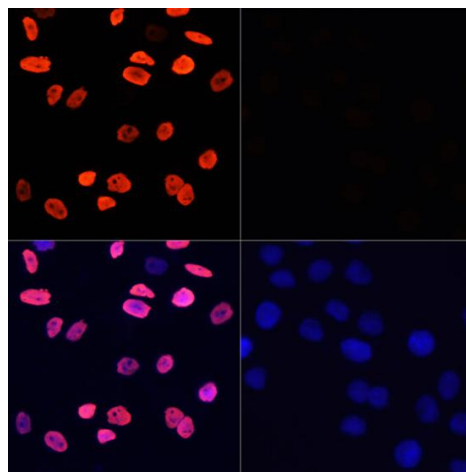
Immunohistochemistry analysis of paraffin-embedded Human colon using Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



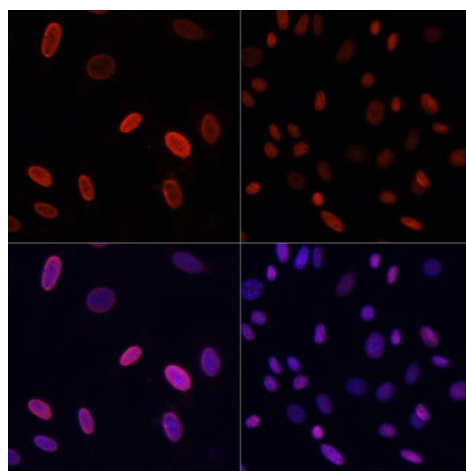
Immunohistochemistry analysis of paraffin-embedded Rat ovary using Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



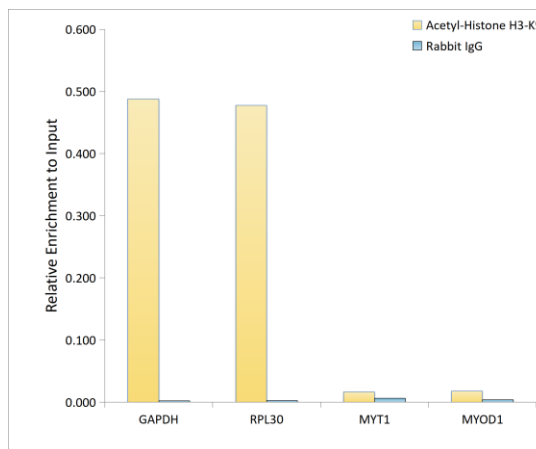
Immunohistochemistry analysis of paraffin-embedded Mouse brain using Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



Immunofluorescence analysis of HeLa cells using Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody at dilution of 1:100 (40x lens). HeLa cells were treated by TSA (1 uM) at 37°C for 18 hours (left). Blue: DAPI for nuclear staining.



Immunofluorescence analysis of NIH/3T3 cells using Histone H3K9ac (Acetyl H3K9) Polyclonal Antibody at dilution of 1:100 (40x lens). NIH/3T3 cells were treated by TSA (1 uM) at 37°C for 18 hours (left). Blue: DAPI for nuclear staining.



Chromatin immunoprecipitation analysis of extracts of HeLa cells, using Histone H3K9ac (Acetyl H3K9) 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.