
Histone H3K9ac (Acetyl H3K9) Monoclonal Antibody [9D2E4G]

(Catalog # A-4055)

Background

Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. This structure consists of approximately 146 bp of DNA wrapped around a nucleosome, an 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 replication-dependent histone that is a member of the histone H3 family. Transcripts from this gene lack polyA tails; instead, they contain a palindromic termination element. This gene is found in the large histone gene cluster on chromosome 6p22-p21.3.

Description

Histone H3K9ac (Acetyl H3K9) Monoclonal Antibody, clone 9D2E4G. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P68431

Purification

Affinity Purified

Immunogen

A synthetic acetylated peptide around K9 of human Histone H3 (NP_003520.1).

Storage

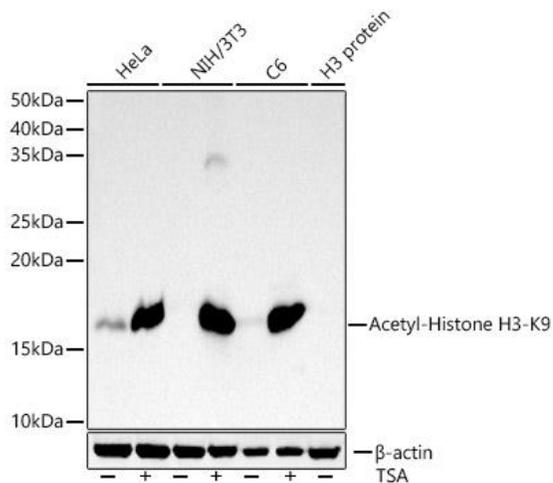
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

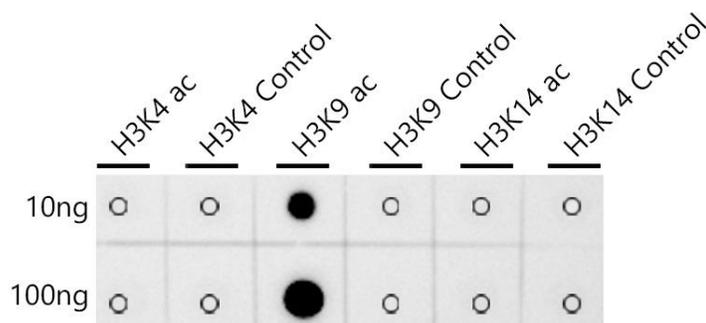
H3/A; H3C2; H3C3; H3C4; H3C6; H3C7; H3C8; H3FA; H3C10; H3C11; H3C12; HIST1H3A

Application

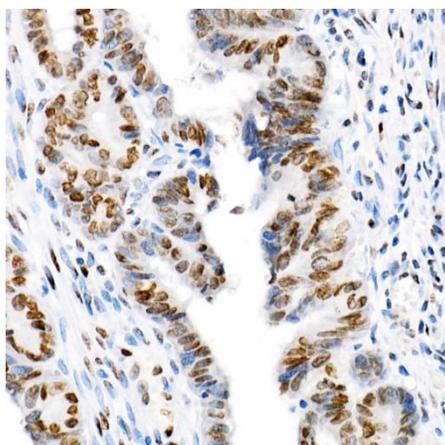
WB, IHC, ChIP, DB; Recommended dilution: DB 1:500 - 1:2000, WB 1:500 - 1:2000, IHC 1:50 - 1:200, ChIP 1:50 - 1:200



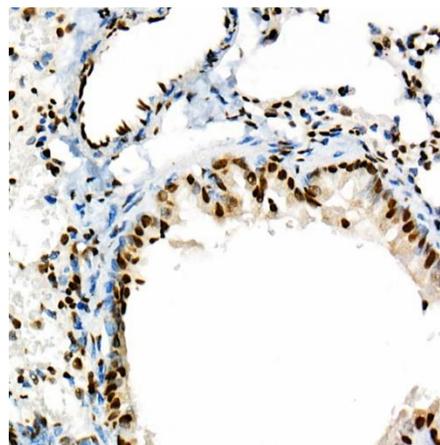
Western blot analysis of extracts of various cell lines, using Histone H3K9ac mAb at 1:1000 dilution. HeLa and NIH3T3, and C6 cells were treated by TSA (1 uM) at 37°C for 18 hours. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25 ug per lane. Blocking buffer: 3% nonfat dry milk in TBST.



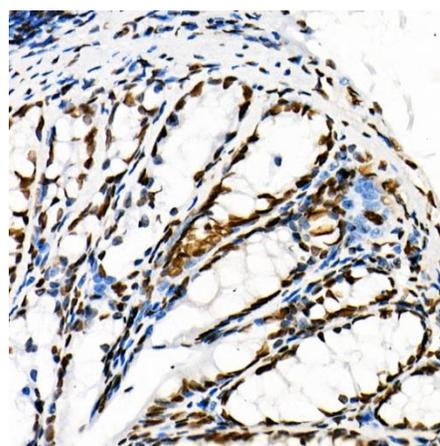
Dot-blot analysis of all sorts of peptides using Histone H3K9ac mAb at 1:1000 dilution.



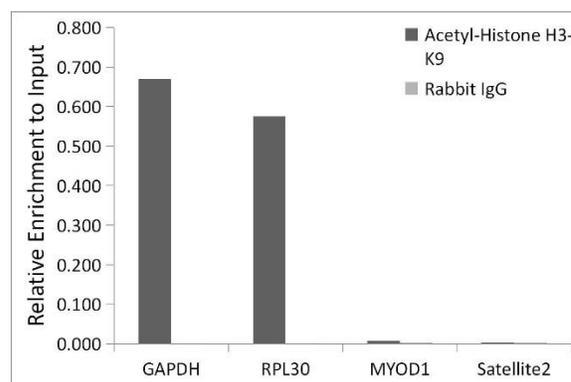
Immunohistochemistry of paraffin-embedded human colon carcinoma using histone Histone H3K9ac mAb at dilution of 1:200 (40x lens).



Immunohistochemistry of paraffin-embedded mouse lung using Histone H3K9ac mAb at dilution of 1:200 (40x lens).



Immunohistochemistry of paraffin-embedded rat colon using Histone H3K9ac mAb at dilution of 1:200 (40x lens).



Chromatin immunoprecipitation analysis of extracts of HeLa cells, using Histone H3K9ac mAb 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.