
Histone H3K14ac (Acetyl H3K14) Polyclonal Antibody

(Catalog # A-4023)

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 H3K14ac (Acetyl H3K14) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.09% Sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat, Wide Range

Isotype

IgG

Uniprot ID

P68431, Q16695

Purification

Affinity Purified

Immunogen

A synthetic acetylated peptide around K14 of human H3 (NP_003520.1).

Storage

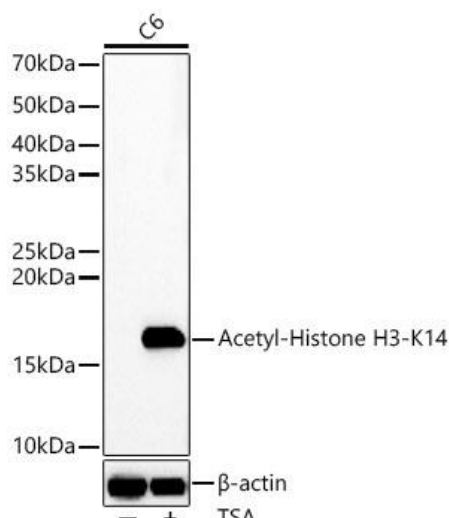
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze/thaw cycles.

Alternative Names

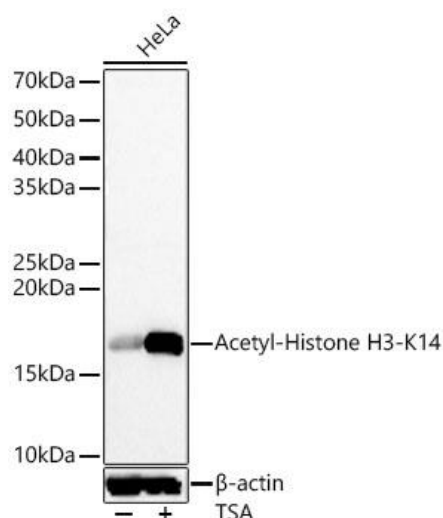
H3K14ac antibody; H3K14a antibody; H3t; H3.4; H3/g; H3FT

Application

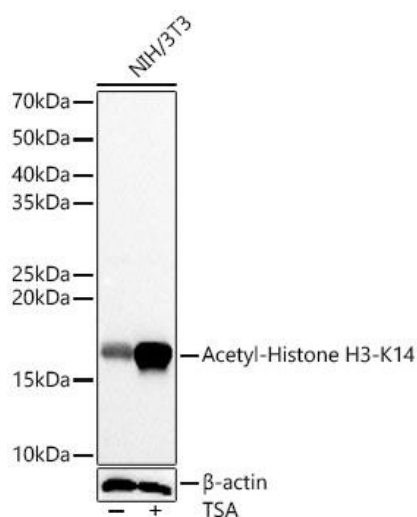
ELISA, WB, IF/ICC; Recommended dilution: WB 1:500-1:1000, IF/ICC 1:50-1:200



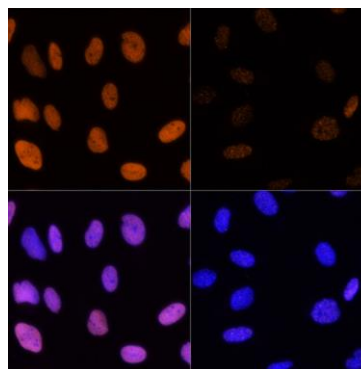
Western blot analysis of lysates from C6 cells using Histone H3K14ac (Acetyl H3K14) Polyclonal Antibody at 1:400 dilution. C6 cells were treated by TSA (1 μ M) at 37°C for 18 hours. Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25 μ g per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 30s.



Western blot analysis of lysates from HeLa cells using Histone H3K14ac (Acetyl H3K14) Polyclonal Antibody at 1:400 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: 30s.



Western blot analysis of lysates from NIH/3T3 cells using Histone H3K14ac (Acetyl H3K14) Polyclonal Antibody at 1:400 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: 30s.



Immunofluorescence analysis of U-2 OS cells using Histone H3K14ac (Acetyl H3K14) Polyclonal Antibody at dilution of 1:100 (40x lens). U-2 OS cells were treated by TSA (1 μ M) at 37°C for 18 hours. Secondary antibody: Cy3-conjugated Goat anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.