
Histone H3K27ac (Acetyl H3K27) Polyclonal Antibody

(Catalog # A-4708)

Histones are proteins found in eukaryotic cell nuclei that package and order the DNA into structural units called nucleosomes. Nucleosomes consist of about 146-147 bp of DNA wrapped around an octamer of histone proteins (histone 2A, histone 2B, histone 3, and histone 4). The N-terminal tails of histones protrude from the globular nucleosome core and can undergo several different types of epigenetic modifications that influence cellular processes. The interaction of a linker histone, H1, with DNA between nucleosomes, facilitates the compaction of chromatin into higher-order structures. This gene is without introns and encodes a histone H3 family member. Transcripts from this gene are missing a polyA tail. As an alternative, they contain palindromic termination elements. This gene is located independently from the other H3 genes. Most H3 genes are found in the histone gene cluster on chromosome 6p22-p21. 3.

Description

Histone H3K27ac (Acetyl H3K27) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Specificity

Broad Range, Human, Mouse, Rat

Formulation

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

Isotype

IgG

Uniprot ID

P68431

Purification

Affinity Purified

Immunogen

A synthetic acetylated peptide around K27 of human Histone H3 (NP_003520.1).

Storage

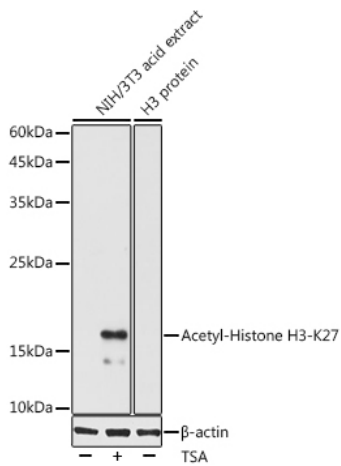
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

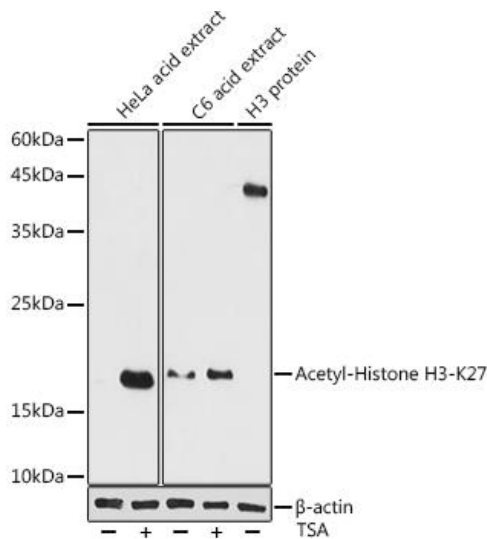
H3t, H3.4, H3/g, H3FT, H3K27ac

Application

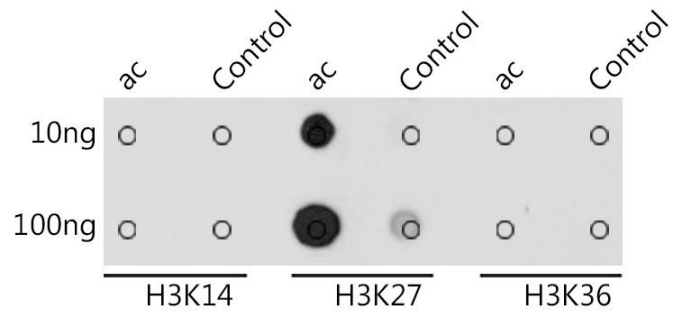
WB, ChIP, ELISA, DB: Recommended dilution, WB 1:500 - 1:1000, ChIP 1:50 - 1:200, DB 1:500 - 1:1000



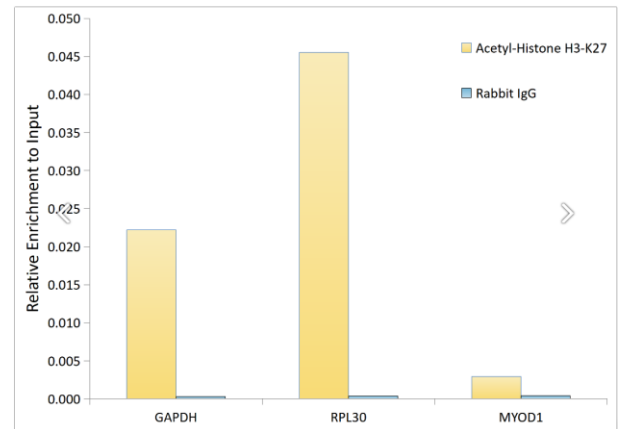
Western blot analysis of extracts of various cell lines, using Acetyl-Histone H3-K27 antibody at 1:1000 dilution. NIH/3T3 cells were treated by TSA (1 μ M) at 37°C for 18 hours. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25 μ g per lane. Blocking buffer: 3% nonfat dry milk in TBST.



Western blot analysis of extracts of various cell lines, using Acetyl-Histone H3-K27 antibody at 1:1000 dilution. HeLa cells and C6 cells were treated by TSA (1 μ M) at 37°C for 18 hours. Secondary antibody: HRP 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.



Western blot analysis of extracts of various cell lines, using Acetyl-Histone H3-K27 antibody at 1:1000 dilution. HeLa cells and C6 cells were treated by TSA (1 μ M) at 37°C for 18 hours. Secondary antibody: HRP 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.



Chromatin immunoprecipitation analysis of extracts of HeLa cells, using Acetyl-Histone H3-K27 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.