
Histone H3K56ac (Acetyl H3K56) Polyclonal Antibody

(Catalog # A-4026)

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, 23, and can also be acetylated at lysine 56.

Description

Histone H3K56ac (Acetyl H3K56) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thimerosal, 50% glycerol, pH7.3.

Specificity

Mouse, Rat, Human, Broad Range

Isotype

IgG

Uniprot ID

Q16695

Purification

Affinity Purified

Immunogen

A synthetic acetylated peptide around K56 of human Histone H3 (NP_003520.1)

Storage

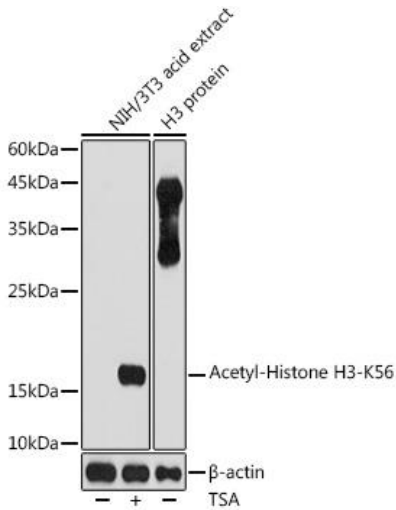
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze.

Alternative Names

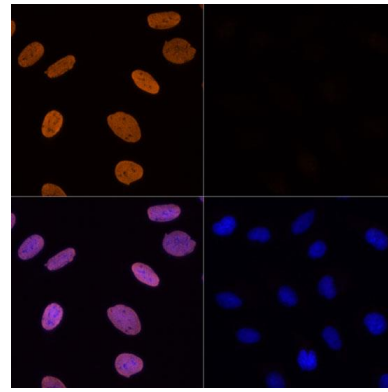
H3K56ac antibody, H3K56a antibody

Application

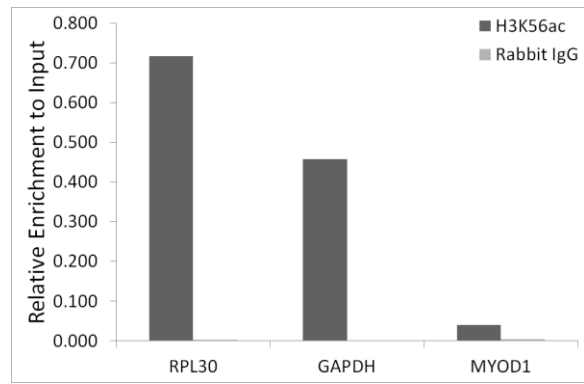
WB, IHC, IF, IP, ChIP, ChIP-seq; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:200, IF 1:50 - 1:200, IP 1:50 - 1:200, ChIP 1:20 - 1:100, ChIP-seq 1:20 - 1:100



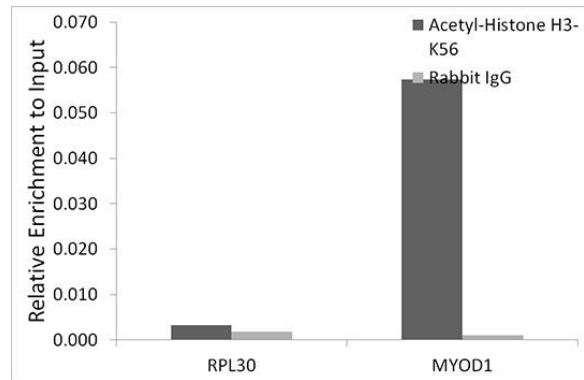
Western blot analysis of extracts of NIH/3T3 cells, using Acetyl-Histone H3-K56 antibody at 1:1000 dilution. NIH/3T3 cells were treated by TSA (1 μ M) at 37°C for 18 hours. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25 μ g per lane. Blocking buffer: 3% nonfat dry milk in TBST.



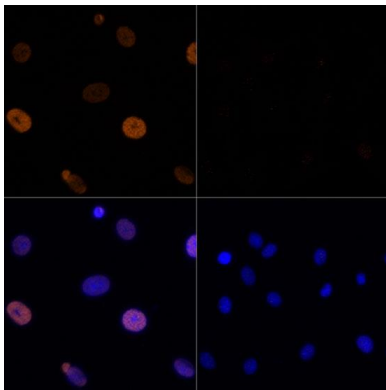
Immunofluorescence analysis of U-2 OS cells using Acetyl-Histone H3-K56 Rabbit pAb at dilution of 100 (40x lens). U-2 OS cells were treated by TSA (1 μ M) at 37°C for 18 hours. Blue: DAPI for nuclear staining.



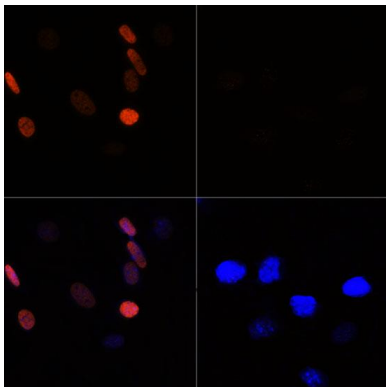
Chromatin immunoprecipitation analysis of extracts of HeLa cells, using Acetyl-Histone H3-K56 antibody and rabbit IgG. The amount of immunoprecipitated DNA was checked by quantitative PCR. Histogram was constructed by the ratios of the immunoprecipitated DNA to the input.



Chromatin immunoprecipitation analysis of extracts of HeLa cells, using Acetyl-Histone H3-K56 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.



Immunofluorescence analysis of C6 cells using Acetyl-Histone H3-K56 Rabbit pAb at dilution of 100 (40x lens). C6 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 Acetyl-Histone H3-K56 Rabbit pAb at dilution of 100 (40x lens). NIH/3T3 cells were treated by TSA (1 μ M) at 37°C for 18 hours. Blue: DAPI for nuclear staining.