

Histone H3K23ac (Acetyl H3K23) Polyclonal Antibody

(Catalog #A-4025)

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 H3K23ac (Acetyl H3K23) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid in PBS containing 50% glycerol, 0.5% BSA* and 0.02% sodium azide.

Specificity

Mouse, Rat, Human

Isotype

IgG

Uniprot ID

P68431/Q71DI3/P84243

Purification

Affinity Purified

Immunogen

Synthesized peptide derived from Human Histone H3 around the acetylation site of K23

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

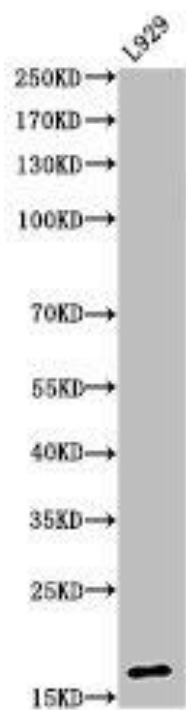
Alternative Names

H3K23ac antibody, H3K23a antibody

Application

WB, IF, ELISA; Recommended dilution: WB:1:500-1:2000, IF:1:200-1:1000, ELISA:1:5000

**The BSA is derived from animal or animal-derived material of negligible amounts. The animal-derived material is subject to heat treatment at a temperature higher than 65°C for at least three hours and acid treatment with pH value of less than 5 for at least three hours, thereby being free of Bovine Spongiform Encephalopathy (BSE) and Transmissible Spongiform Encephalopathy (TSE). Additionally, as all BSA-containing products are strictly intended for research purposes and not for diagnostic or therapeutic use, they are not subject to certification authority oversight.*



Western Blot analysis of L929 cells using Histone H3K23ac (Acetyl H3K23) Polyclonal Antibody.



Western Blot analysis of L929 cells using Histone H3K23ac (Acetyl H3K23) Polyclonal Antibody.