

---

## 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. Aliquot store at -20°C . Avoid repeated freeze / thaw cycles.

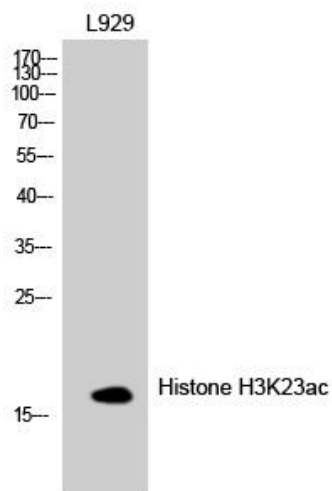
### Alternative Names

H3K23ac antibody, H3K23a antibody

### Application

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

---



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