

---

## Histone H4-K16ac (Acetyl H4K16) Polyclonal Antibody

(Catalog # A68398)

---

### Background

Histone H4- along with H2A, H2B, and H3- is involved in the structure of chromatin in eukaryotic cells. Histone H4 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 H4, which results in remodeling of the nucleosome structure into an open conformation more accessible to transcription complexes. In most species, histone H4 is primarily acetylated at lysine 5, 8, 12, and 16.

### Description

Histone H4-K16ac (Acetyl H4K16) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

P62805

### Purification

Affinity Purified

### Immunogen

Synthesized peptide derived from Human Histone H4 around the acetylation site of K16

### Storage

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

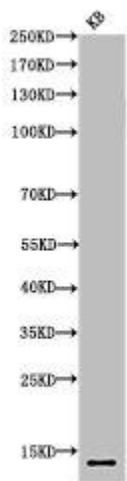
### Alternative Names

HIST1H4A; H4/A; H4FA; HIST1H4B; H4/I; H4FI; HIST1H4C; H4/G; H4FG; HIST1H4D; H4/B; H4FB; HIST1H4E; H4/J; H4FJ; HIST1H4F; H4/C; H4FC; HIST1H4H; H4/H; H4FH; HIST1H4I; H4/M; H4FM; HIST1H4J; H4/E; H4FE; HIST1H4K; H4/D; H4FD; HIST1H4L; H4/K; H4FK

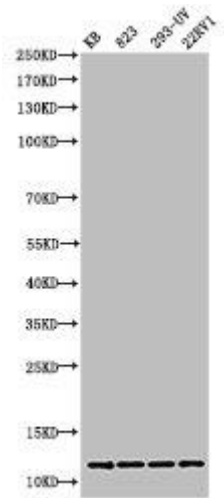
### Application

WB, IHC, IF, ELISA; Recommended dilution: WB:1:500-1:2000, IHC:1:100-1:300, IF:1:200-1:1000, ELISA:1:5000

---



Western Blot analysis of KB cells using Histone H4-K16ac (Acetyl H4K16) Polyclonal Antibody



Western Blot analysis of KB 823 293-UV 22RV1 cells using Histone H4-K16ac (Acetyl H4K16) Polyclonal Antibody