

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

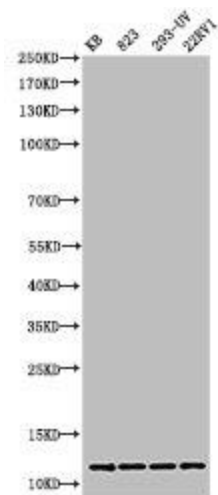
Application

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

**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 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