

Histone H4K12ac (Acetyl H4K12) Polyclonal Antibody

(Catalog #A-4029)

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 H4K12ac (Acetyl H4K12) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Purification

Affinity Purified

Immunogen

Synthesized peptide derived from human Histone H4 around the acetylation site of K12.

Specificity

Human, Mouse, Rat, Monkey

Uniprot No

P62805

Isotype

IgG

Formulation

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

Storage

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

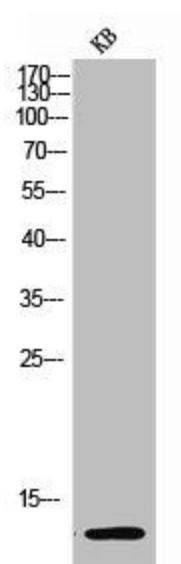
Alternative Names

H4K12ac antibody, H4K12a antibody

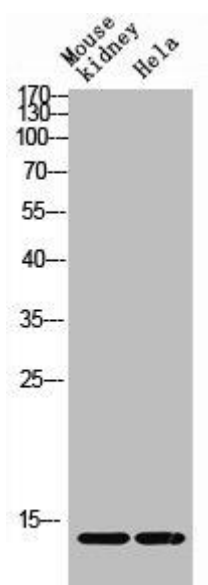
Applications

ELISA, IF, IHC, WB; Recommended dilution: WB:1:500-1:2000, IHC:1:100-1:300, IF:1:200-1:1000, ELISA:1:10000

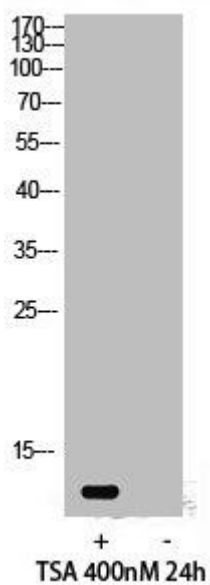
**The BSA is derived from animal or animal-derived material of negligible amounts. The animal-derived material is subject to heat treatment at a temperature higher than 65°C for at least three hours and acid treatment with pH value of less than 5 for at least three hours, thereby being free of Bovine 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 H4K12ac (Acetyl H4K12) Polyclonal Antibody.



Western blot analysis of Mouse-kidney HELA lysis using Histone H4K12ac (Acetyl H4K12) Polyclonal Antibody.



Western Blot analysis of 293 cells using Histone H4K12ac (Acetyl H4K12) Polyclonal Antibody.