
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

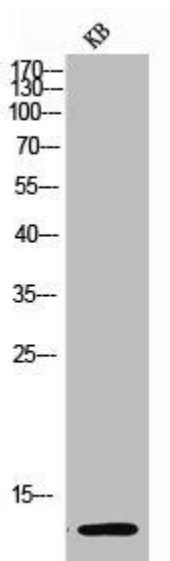
Store at -20°C or -80°C. Avoid repeated freeze

Alternative Names

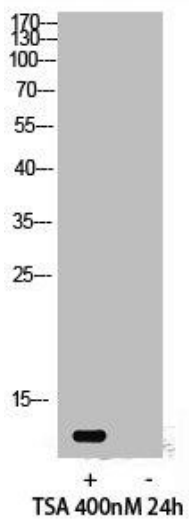
H4K12ac antibody, H4K12a antibody

Application

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



Western Blot analysis of KB cells using Acetyl-Histone H4 (K12) Polyclonal Antibody



Western Blot analysis of 293 cells using Acetyl-Histone H4 (K12) Polyclonal Antibody