

Histone H4K8ac (Acetyl H4K8) Polyclonal Antibody

(Catalog #A-4028)

Background

Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. This structure consists of approximately 146 bp of DNA wrapped around a nucleosome, an octamer composed of pairs of each of the four core histones (H2A, H2B, H3, and H4). The chromatin fiber is further compacted through the interaction of a linker histone, H1, with the DNA between the nucleosomes to form higher order chromatin structures. This gene is intronless and encodes a member of the histone H4 family. Transcripts from this gene lack polyA tails; instead, they contain a palindromic termination element. This gene is found in a histone cluster on chromosome 1. This gene is one of four histone genes in the cluster that are duplicated; this record represents the centromeric copy.

Description

Histone H4K8ac (Acetyl H4K8) Polyclonal Antibody. Unconjugated. Raised in: Rabbit

Purification

Affinity purified

Specificity

Human, Mouse, Rat, Monkey

Isotype

IgG

Immunogen

Synthesized peptide derived from human Histone H4 around the acetylation site of K8.

Formulation

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

Storage

Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

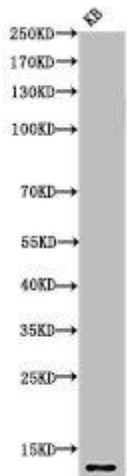
Alternative Names

H4K8ac antibody; H4K8a antibody; H4; H4/n; H4F2; H4FN; FO108; HIST2H4

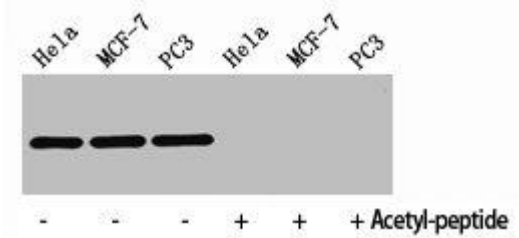
Application

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 H4K8ac (Acetyl H4K8) Polyclonal Antibody



Western Blot analysis of HELA MCF7 PC3 cells using Histone H4K8ac (Acetyl H4K8) Polyclonal Antibody