

Histone H4K12ac (Acetyl H4K12) Polyclonal Antibody

(Catalog # A68357)

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.

Formulation

PBS with 0.02% sodium azide, 50% glycerol, pH 7.4.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P62805

Immunogen

Synthetic Peptide

Storage

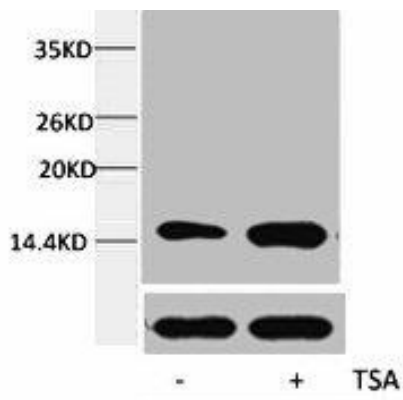
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

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; HIST2H4A; H4/

Application

WB: 1:1000-2000



Western blot analysis of extracts from Hela cells, untreated (-) or treated, at a 1:5000 dilution using the Histone H4K12ac (Acetyl H4K12) Polyclonal Antibody.