
Histone H3K14ac (Acetyl H3K14) Polyclonal Antibody

(Catalog # A-4023)

Background

Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. Nucleosomes consist of approximately 146 bp of DNA wrapped around a histone 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 H3 family. Transcripts from this gene lack polyA tails; instead, they contain a palindromic termination element. This gene is located separately from the other H3 genes that are in the histone gene cluster on chromosome 6p22-p21. 3.

Description

Histone H3K14ac (Acetyl H3K14) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P68431, Q71DI3, P84243

Purification

Affinity Purified

Immunogen

Synthesized peptide derived from Human Histone H3 around the acetylation site of K14.

Storage

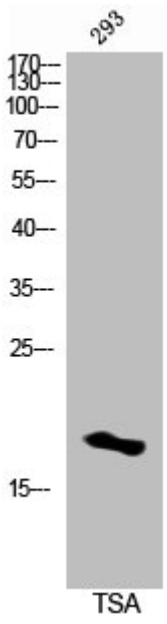
Shipped at 4°C. Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

H3K14ac antibody; H3K14a antibody; H3t; H3.4; H3/g; H3FT

Application

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



Western Blot analysis of 3T3 cells using Histone H3K14ac Polyclonal Antibody