
Histone H4K20ac (Acetyl H4K20) Polyclonal Antibody

(Catalog # A73223)

Background

Core component of nucleosome. Nucleosomes wrap and compact DNA into chromatin, limiting DNA accessibility to the cellular machineries which require DNA as a template. Histones thereby play a central role in transcription regulation, DNA repair, DNA replication and chromosomal stability. DNA accessibility is regulated via a complex set of post-translational modifications of histones, also called histone code, and nucleosome remodeling.

Description

Histone H4K20ac (Acetyl H4K20) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300; 50% Glycerol, 0.01M PBS, pH 7.4

Specificity

Human

Isotype

IgG

Uniprot ID

P62805

Purification

Affinity Purified

Immunogen

Peptide sequence around site of Acetyl-Lys (20) derived from Human Histone H4

Storage

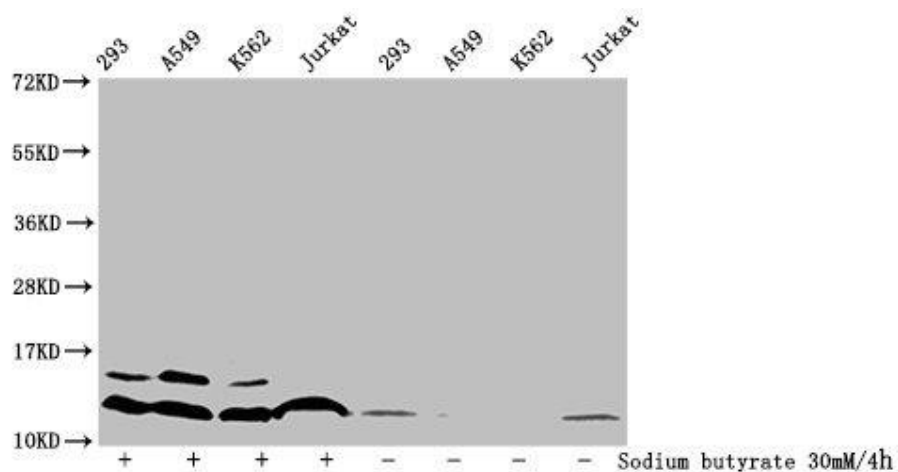
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Acetyl-HIST1H4A (K20); Histone H4, HIST1H4A; HIST1H4B; HIST1H4C; HIST1H4D; HIST1H4E; HIST1H4F; HIST1H4H; HIST1H4I; HIST1H4J; HIST1H4K; HIST1H4L; HIST2H4A; HIST2H4B; HIST4H4, H4/A H4FA; H4/I H4FI; H4/G H4FG; H4/B H4FB; H4/J H4FJ; H4/C H4FC; H4/H H4FH; H4/M H4FM; H4/E H4FE; H4/D H4FD; H4/K H4FK; H4/N H4F2 H4FN HIST2H4; H4/O H4FO;

Application

ELISA, WB, IF; Recommended dilution: WB:1:100-1:1000, IF:1:1-1:10



Western Blot

Detected samples: 293 whole cell lysate, A549 whole cell lysate, K562 whole cell lysate, Jurkat whole cell lysate; Untreated (-) or treated (+) with 30mM sodium butyrate for 4h

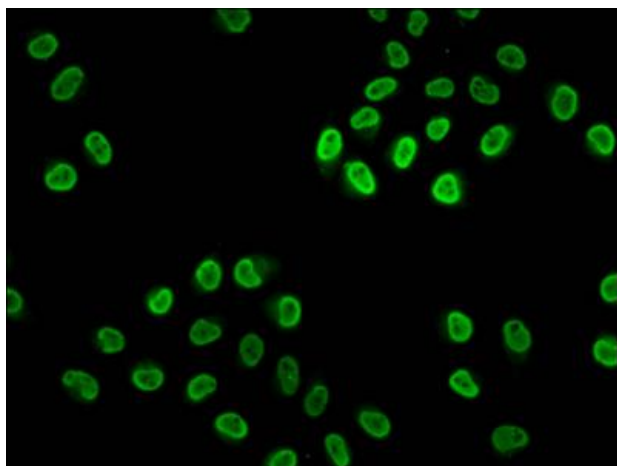
All lanes: Histone H4K20ac (Acetyl H4K20) Polyclonal Antibody at 1:100

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 12 kDa

Observed band size: 12 kDa



Immunofluorescence staining of HeLa cells (treated with 30mM sodium butyrate for 4h) with Histone H4K20ac (Acetyl H4K20) Polyclonal Antibody at 1:2.5, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C.