

Histone H4K79ac (Acetyl H4K79) Polyclonal Antibody

(Catalog # A-4053)

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 H4K79ac (Acetyl H4K79) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. Preservative: 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 (79) derived from Human Histone H4

Storage

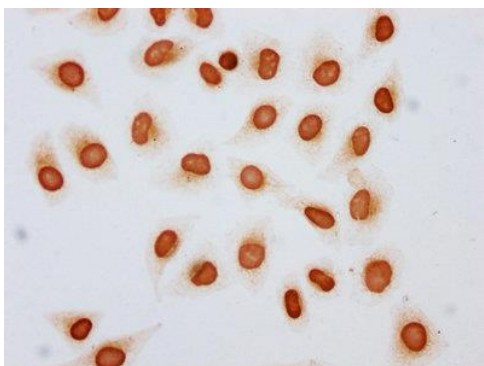
Shipped at 4°C (Blue Ice). Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

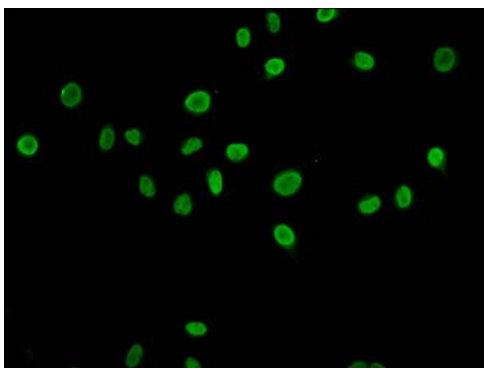
H4 antibody; H4.k antibody; H4/a antibody; H4/b antibody; H4/c antibody; H4/d antibody; H4/e antibody; H4/g antibody; H4/h antibody; H4/l antibody; H4/j antibody; H4/k antibody; H4/m antibody; H4/n antibody; H4/p antibody; H4_HUMAN antibody; H4F2 antibody; H4F2iii antibody; H4F2iv antibody; H4FA antibody; H4FB antibody; H4FC antibody; H4FD antibody; H4FE antibody; H4FG antibody; H4FH antibody; H4FI antibody; H4FJ antibody; H4FK antibody; H4FM antibody; H4FN antibody; H4M antibody; HIST1H4A antibody; HIST1H4B antibody; HIST1H4C antibody; HIST1H4D antibody; HIST1H4E antibody; HIST1H4F antibody; HIST1H4H antibody; HIST1H4I antibody; HIST1H4J antibody; HIST1H4K antibody; HIST1H4L antibody; HIST2H4 antibody; HIST2H4A antibody; Hist4h4 antibody; Histone 1 H4a antibody; Histone 1 H4b antibody; Histone 1 H4c antibody; Histone 1 H4d antibody; Histone 1 H4e antibody; Histone 1 H4f antibody; Histone 1 H4h antibody; Histone 1 H4i antibody; Histone 1 H4k antibody; Histone 1 H4l antibody; Histone 2 H4a antibody; histone H4 antibody;

Application

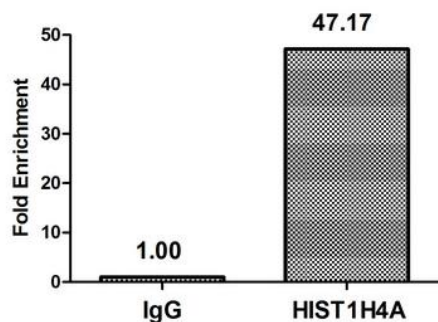
ELISA, ICC, IF, ChIP; Recommended dilution: ICC 1:10-1:100, IF 1:1-1:10



Immunocytochemistry analysis of Histone H4K79ac (Acetyl H4K79) Polyclonal Antibody diluted at 1:10 and staining in HeLa cells (treated with 30mM sodium butyrate for 4h) was performed on a Leica Bond™ system. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Immunofluorescence staining of HeLa cells (treated with 30mM sodium butyrate for 4h) with Histone H4K79ac (Acetyl H4K79) Polyclonal Antibody at 1:5, 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. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).



Chromatin Immunoprecipitation HeLa (4×10^6 , treated with 30mM sodium butyrate for 4h) were treated with Benzanase, sonicated, and immunoprecipitated with 5µg of Histone H4K79ac (Acetyl H4K79) Polyclonal Antibody or a control normal rabbit IgG. The resulting ChIP DNA was quantified using real-time PCR with primers against the beta-Globin promoter.