

Histone H4 Acetylation Antibody Panel Pack

Base Catalog # C10013

PACK CONTENTS

Component	Size	Shipping Temperature	Storage Upon Receipt	Storage Checklist
4K5A Histone H4K5ac (Acetyl H4K5) Polyclonal Antibody	25 µl	4°C	–20°C	
4K8A Histone H4K8ac (Acetyl H4K8) Polyclonal Antibody	25 µl	4°C	–20°C	
4K12A Histone H4K12ac (Acetyl H4K12) Polyclonal Antibody	25 µl	4°C	–20°C	
4K16A Histone H4K16ac (Acetyl H4K16) Polyclonal Antibody	25 µl	4°C	–20°C	
HGR2 HRP-Goat Anti-Rabbit Secondary Antibody	50 µg	4°C	–20°C	

SHIPPING & STORAGE

This product is shipped on frozen ice packs at 4°C. Upon receipt: (1) Store all components at –20°C away from light.

All components of the product are stable for 6 months from the date of shipment, when stored properly.

Histone H4K5ac (Acetyl H4K5) Polyclonal Antibody

Component Cat. #C10013-1-4K5A

Description

Histone H4K5ac (Acetyl H4K5) 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

Purification

Affinity Purified

Immunogen

Synthetic Peptide

Storage

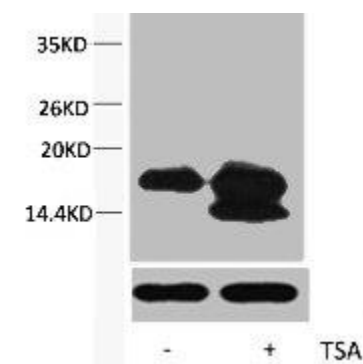
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

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; Recommended dilution: WB: 1:2000-5000



Western blot analysis of extracts from Hela cells, untreated (-) or treated, 1:5000.

Histone H4K8ac (Acetyl H4K8) Polyclonal Antibody

Component Cat. #C10013-1-4K8A

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 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

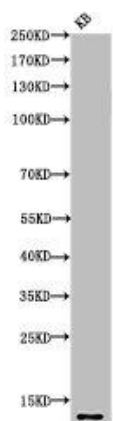
Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

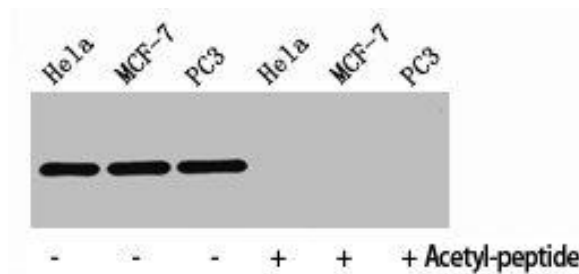
H4K8ac antibody; H4K8a antibody; H4; H4/n; H4F2; H4FN; FO108; HIST2H4

Application

WB, IHC, IF, ELISA, ChIP; Recommended dilution: WB:1/500-1/2000, IHC:1/100-1/300, IF:1/200-1/1000, ELISA:1/5000



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.

Histone H4K12ac (Acetyl H4K12) Polyclonal Antibody

Component Cat. #C10013-1-4K12A

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.

Purification

Affinity Purified

Immunogen

Synthesized peptide derived from human Histone H4 around the acetylation site of K12.

Specificity

Human, Mouse, Rat, Monkey

Uniprot No

P62805

Isotype

IgG

Formulation

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

Storage

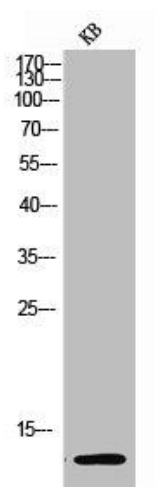
Store at -20°C or -80°C. Avoid repeated freeze

Alternative Names

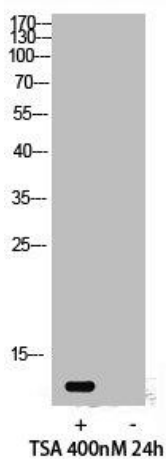
H4K12ac antibody, H4K12a antibody

Application

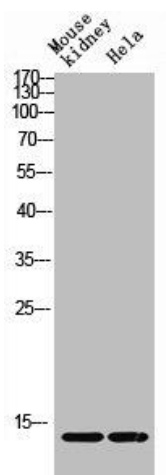
ELISA, IF, IHC, WB; Recommended dilution: WB:1:500-1:2000, IHC:1:100-1:300, IF:1:200-1:1000, ELISA:1:10000



Western Blot analysis of KB cells using Acetyl-Histone H4 (K12) Polyclonal Antibody



Western Blot analysis of 293 cells using Acetyl-Histone H4 (K12) Polyclonal Antibody



Western blot analysis of Mouse-kidney HELA lysis using Acetyl-Histone H4 (K12) antibody.

Histone H4-K16ac (Acetyl H4K16) Polyclonal Antibody

Component Cat. #C10013-1-4K16A

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 H4-K16ac (Acetyl H4K16) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P62805

Purification

Affinity Purified

Immunogen

Synthesized peptide derived from Human Histone H4 around the acetylation site of K16

Storage

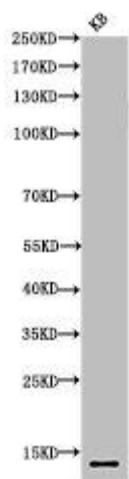
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

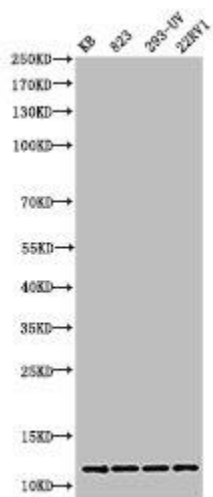
HIST1H4A; H4/A; H4FA; HIST1H4B; H4/I; H4FI; HIST1H4C; H4/G; H4FG; HIST1H4D; H4/B; H4FB; HIST1H4E; H4/J; H4FJ; HIST1H4F; H4/C; H4FC; HIST1H4H; H4/H; H4FH; HIST1H4I; H4/M; H4FM; HIST1H4J; H4/E; H4FE; HIST1H4K; H4/D; H4FD; HIST1H4L; H4/K; H4FK

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



Western Blot analysis of KB cells using Histone H4-K16ac (Acetyl H4K16) Polyclonal Antibody



Western Blot analysis of KB 823 293-UV 22RV1 cells using Histone H4-K16ac (Acetyl H4K16) Polyclonal Antibody

HRP- Goat Anti-Rabbit Secondary Antibody

Component Cat. #C10013-1-HGR2

Description

Goat anti-rabbit IgG recognizes rabbit IgG whole molecule. This secondary antibody was purified using antigen affinity chromatography. The antibody is conjugated with peroxidase.

Antibody Type

Polyclonal Antibody

Purification

Liquid; this product was prepared from monospecific antiserum by immunoaffinity chromatography, followed by solid phase adsorption(s) to remove any unwanted reactivities.

Immunogen

Rabbit IgG whole molecule

Isotype

IgG

Formulation

In 10 mM sodium phosphate, 75 mM NaCl, 50% (v/v) glycerol, pH 7.2.

Specificity

Rabbit

Storage

Store at -20°C. Aliquot to avoid repeated freezing and thawing.

Handling Recommendations

The optimal working dilution should be determined by the end user. For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions

Western Blot: 1:1000-1: 10000; Immunohistochemistry: 1:100-1:500; Immunofluorescence: 1:100-1:500; ELISA: 1:2000-1:20000

RELATED PRODUCTS

Histone Modification Antibodies

A68356	Histone H4K5ac (Acetyl H4K5) Polyclonal Antibody
A-4028	Histone H4K8ac (Acetyl H4K8) Polyclonal Antibody
A-4029	Histone H4K12ac (Acetyl H4K12) Polyclonal Antibody
A68398	Histone H4K16ac (Acetyl H4K16) Polyclonal Antibody
A12004	HRP-Goat Anti-Rabbit Secondary Antibody

Histone Modification Panel Packs

C10000	Histone H3 Methylation Antibody Panel Pack I – Active Genes
C10001	Histone H3 Methylation Antibody Panel Pack I – Repression Genes
C10002	Histone H3 Methylation Antibody Panel Pack II – Active Genes
C10003	Histone H3 Methylation Antibody Panel Pack II – Repression Genes
C10004	Histone H3 Methylation Antibody Panel Pack III – Active Genes
C10005	Histone H3K4 Methylation Antibody Panel Pack
C10006	Histone H3K9 Methylation Antibody Panel Pack
C10007	Histone H3K27 Methylation Antibody Panel Pack
C10008	Histone H3K36 Methylation Antibody Panel Pack
C10009	Histone H3K79 Methylation Antibody Panel Pack
C10010	Histone H3 Acetylation Antibody Panel Pack I
C10011	Histone H3 Acetylation Antibody Panel Pack II
C10012	Histone H4K20 Methylation Antibody Panel Pack
C10013	Histone H4 Acetylation Antibody Panel Pack
C10014	Histone H3 Phosphorylation Antibody Panel Pack
C10015	Histone H3R2 Methylation Antibody Panel Pack
C10016	Histone H3R8 Methylation Antibody Panel Pack
C10017	Histone H3R17 Methylation Antibody Panel Pack
C10018	Histone H3R26 Methylation Antibody Panel Pack
C10019	Histone H4R3 Methylation Antibody Panel Pack