

# Histone H3 Acetylation Antibody Panel Pack II

Base Catalog # C10011

## PACK CONTENTS

Component	Size	Shipping Temperature	Storage Upon Receipt	Storage Checklist
<b>3K23A</b> Histone H3K23ac (Acetyl H3K23) Polyclonal Antibody	25 µl	4°C	-20°C	
<b>3K36A</b> Histone H3K36ac (Acetyl H3K36) Polyclonal Antibody	25 µl	4°C	-20°C	
<b>3K56A</b> Histone H3K56ac (Acetyl H3K56) Polyclonal Antibody	25 µl	4°C	-20°C	
<b>3K79A</b> Histone H3K79ac (Acetyl H3K79) Polyclonal Antibody	25 µl	4°C	-20°C	
<b>HGR2</b> 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) 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 H3K23ac (Acetyl H3K23) Polyclonal Antibody

Component Cat. #C10011-3K23A

### Background

Histone H3 along with H2A, H2B and H4 is involved in the structure of chromatin in eukaryotic cells. Histone H3 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 H3, which results in remodeling of the nucleosome structure into an open conformation more accessible to transcription complexes. In most species, histone H3 is primarily acetylated at lysine 9, 14, 18, and 23.

### Description

Histone H3K23ac (Acetyl H3K23) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Mouse, Rat, Human

### Isotype

IgG

### Uniprot ID

P68431/Q71DI3/P84243

### Purification

Affinity Purified

### Immunogen

Synthesized peptide derived from Human Histone H3 around the acetylation site of K23

### Storage

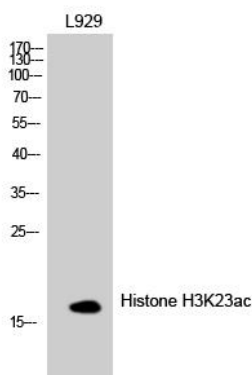
Shipped at 4°C. Aliquot store at -20°C. Avoid repeated freeze / thaw cycles.

### Alternative Names

H3K23ac antibody, H3K23a antibody

### Application

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



Western Blot analysis of L929 cells using Histone H3K23ac (Acetyl H3K23) Polyclonal Antibody.

## Histone H3K36ac (Acetyl H3K36) Polyclonal Antibody

Component Cat. #C10011-3K36A

### Description

Histone H3K36ac (Acetyl H3K36) 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

P68431/Q71DI3/P84243

### Purification

Affinity Purified

### Immunogen

Synthetic Peptide

### Storage

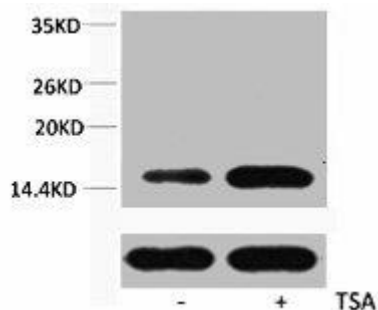
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

HIST1H3A; H3FA; HIST1H3B; H3FL; HIST1H3C; H3FC; HIST1H3D; H3FB; HIST1H3E; H3FD; HIST1H3F; H3FI; HIST1H3G; H3FH; HIST1H3H; H3FK; HIST1H3I; H3FF; HIST1H3J; H3FJ; Histone H3.1; Histone H3/a; Histone H3/b; Histone H3/c; Histone H3/d; Histone H3/f; Histone H3/

### Application

WB; Recommended dilution: WB: 1:200-500



Western blot analysis of extracts from HeLa cells, untreated (-) or treated, 1:2000.

## Histone H3K56ac (Acetyl H3K56) Polyclonal Antibody

Component Cat. #C10011-3K56A

### Background

Histone H3 along with H2A, H2B and H4 is involved in the structure of chromatin in eukaryotic cells. Histone H3 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 H3, which results in remodeling of the nucleosome structure into an open conformation more accessible to transcription complexes. In most species, histone H3 is primarily acetylated at lysine 9, 14, 18, 23, and can also be acetylated at lysine 56.

### Description

Histone H3K56ac (Acetyl H3K56) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Liquid. Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3 .

### Specificity

Mouse, Rat, Human, Broad Range

### Isotype

IgG

### Uniprot ID

Q16695

### Purification

Affinity Purified

### Immunogen

Synthetic Peptide of Human Acetyl-Histone H3-K56

### Storage

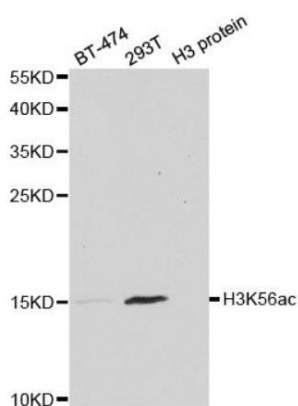
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

### Alternative Names

H3K56ac antibody, H3K56a antibody

### Application

WB, IHC, IF, IP, ChIP, ChIPseq



WB analysis of Histone H3K56ac (Acetyl H3K56) Polyclonal Antibody.

## Histone H3K79ac (Acetyl H3K79) Polyclonal Antibody

Component Cat. #C10011-3K79A

### 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 H3K79ac (Acetyl H3K79) 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

P68431/Q71DI3/P84243

### Purification

Affinity Purified

### Immunogen

Synthetic Peptide

### Storage

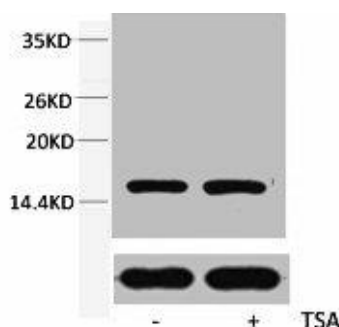
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

HIST1H3A; H3FA; HIST1H3B; H3FL; HIST1H3C; H3FC; HIST1H3D; H3FB; HIST1H3E; H3FD; HIST1H3F; H3FI; HIST1H3G; H3FH; HIST1H3H; H3FK; HIST1H3I; H3FF; HIST1H3J; H3FJ; Histone H3.1; Histone H3/a; Histone H3/b; Histone H3/c; Histone H3/d; Histone H3/f; Histone H3/

### Application

WB, ELISA; Recommended dilution: WB:1:500-1:1000, ELISA:1:20000



Western blot analysis of extracts from HeLa cells, untreated (-) or treated, 1:2000.

---

## HRP- Goat Anti-Rabbit Secondary Antibody

Component Cat. #C10011- 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

A-4025	Histone H3K23ac (Acetyl H3K23) Polyclonal Antibody
A68348	Histone H3K36ac (Acetyl H3K36) Polyclonal Antibody
A-4026	Histone H3K56ac (Acetyl H3K56) Polyclonal Antibody
A68349	Histone H3K79ac (Acetyl H3K79) 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