

Histone H3 Phosphorylation Antibody Panel Pack

Base Catalog # C10014

PACK CONTENTS

| Component | Size | Shipping Temperature | Storage Upon Receipt | Storage Checklist |
|---|-------|----------------------|----------------------|-------------------|
| 3SR10 Phospho Histone H3 (Ser10) Polyclonal Antibody | 25 µl | 4°C | –20°C | |
| 3SR28 Phospho Histone H3 (Ser28) Polyclonal Antibody | 25 µl | 4°C | –20°C | |
| 3TH11 Phospho Histone H3 (Thr11) 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.

Phospho Histone H3 (Ser10) Polyclonal Antibody

Component Cat. #C10014-1-3SR10

Description

Phospho Histone H3 (Ser10) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 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

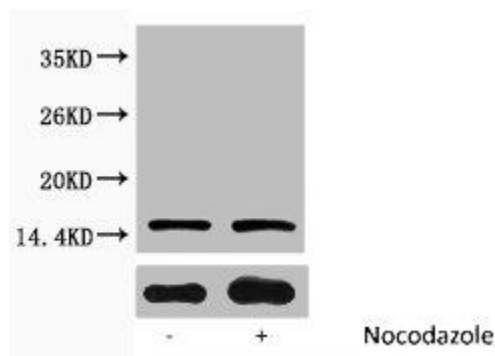
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

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/; H3Ser10; H3Ser10ph; H3S10ph

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.

Phospho-Histone H3 (Ser28) Polyclonal Antibody

Component Cat. #C10014-1-3SR28

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 phosphorylated at serine 10, serine 28, threonine 3 and threonine 11.

Description

Phospho-Histone H3 (Ser28) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. PBS, pH 7.4, containing 0.02% sodium azide as Preservative and 50% Glycerol.

Specificity

Mouse, Rat, Human

Isotype

IgG

Uniprot ID

P68431/Q71DI3/P84243

Purification

Affinity Purified

Immunogen

Synthetic Peptide

Storage

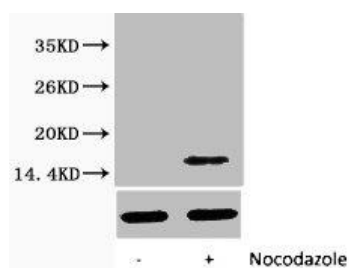
Shipped at 4°C (Blue Ice). Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

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.

Phospho-Histone H3 (Thr11) Polyclonal Antibody

Component Cat. #C10014-1-3TH11

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 phosphorylated at serine 10, serine 28, threonine 3, and threonine 11.

Description

Phospho-Histone H3 (Thr11) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. PBS, pH 7.4, containing 0.02% sodium azide as Preservative and 50% Glycerol.

Specificity

Mouse, Rat, Human

Isotype

IgG

Uniprot ID

P68431/Q71DI3/P84243

Purification

Affinity Purified

Immunogen

Synthetic Peptide

Storage

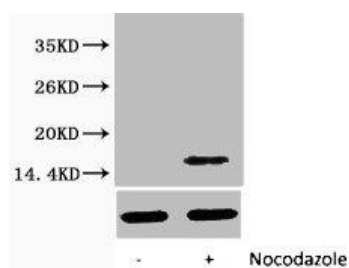
Shipped at 4°C (Blue Ice). Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

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. #C10014-1-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

| | |
|--------|--|
| A68397 | Phospho Histone H3 (Ser10) Polyclonal Antibody |
| A-4050 | Phospho Histone H3 (Ser28) Polyclonal Antibody |
| A-4052 | Phospho Histone H3 (Thr11) 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 |