

## Succinyl HIST1H4A (K12) Polyclonal Antibody

(Catalog # A70383)

### 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

Succinyl HIST1H4A (K12) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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 Succinyl-Lys (12) derived from Human Histone H4

### Storage

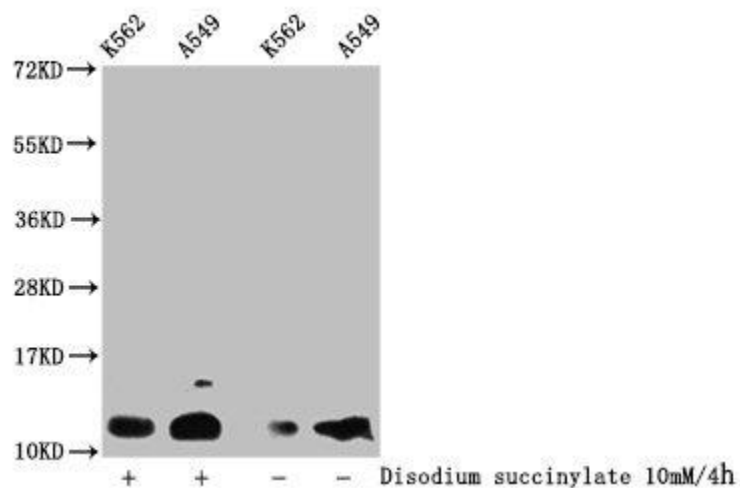
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

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

### Application

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



#### Western Blot

Detected samples: K562 whole cell lysate, A549 whole cell lysate; Untreated (-) or treated (+) with 10mM disodium succinylate for 4h

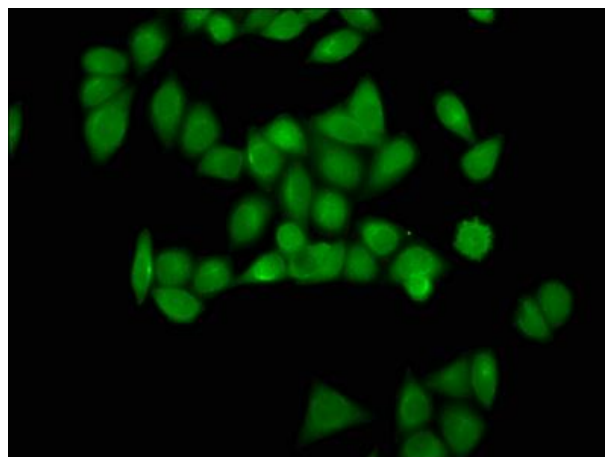
All lanes: Succinyl HIST1H4A (K12) Polyclonal Antibody at 1:100

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 10mM disodium succinylate for 4h) with Succinyl HIST1H4A (K12) Polyclonal Antibody, 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.