

## Phospho Histone H3 (Ser10) Polyclonal Antibody

(Catalog # A68397)

### 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

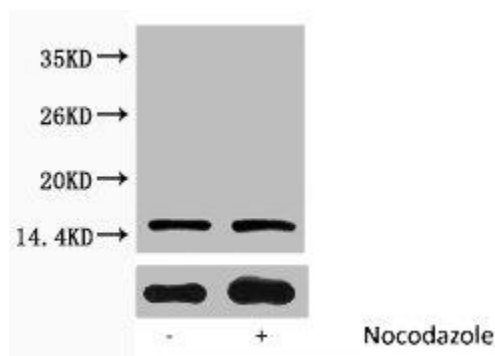
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### 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.