

---

## Histone H3 Monoclonal Antibody [1G1]

(Catalog # A68386)

---

### Description

Histone H3 Monoclonal Antibody [1G1]. Unconjugated. Raised in: Mouse.

### Formulation

PBS with 0.02% sodium azide, 50% glycerol, pH 7.4.

### Specificity

Human, Mouse, Rat, Yeast

### Isotype

IgG

### Uniprot ID

P68431/Q71DI3/P84243

### Purification

Affinity Purified

### Immunogen

Recombinant Protein

### Storage

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

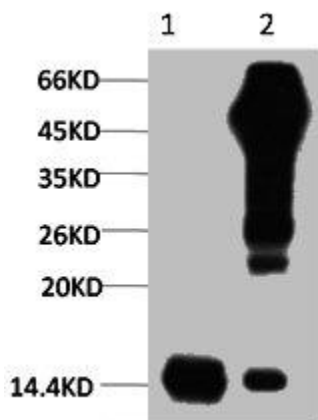
### Alternative Names

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/

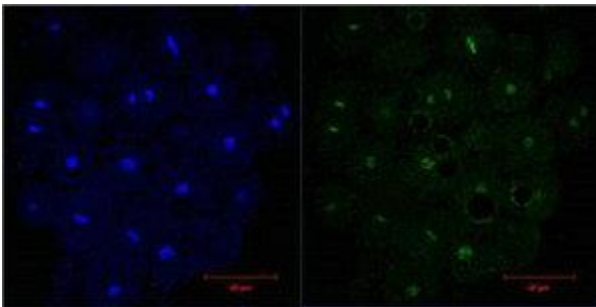
### Application

ELISA, WB, IF, IP; Recommended dilution: WB:1:500-1:5000, IF:1:50-1:200, IP:1:200-1:2000

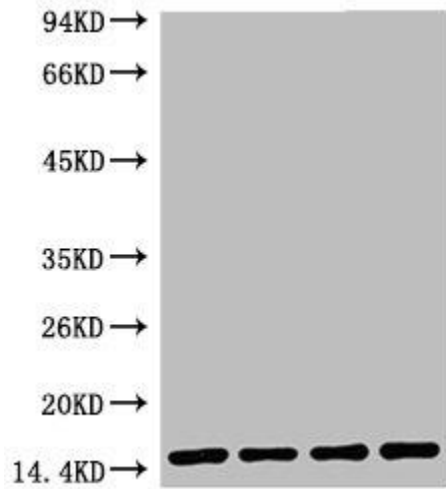
---



1) Input: HeLa Cell Lysate 2) IP product: IP dilute 1:200



IF analysis of HeLa, diluted at 1:200.



Western blot analysis of 1) HeLa, 2) Raw, 3) Mouse Brain Tissue, 4) Rat Brain Tissue, diluted at 1:5000.