

## Histone H4K59me2 (H4K59 Dimethyl) Polyclonal Antibody

(Catalog # A68337)

### Description

Histone H4K59me2 (H4K59 Dimethyl) 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

P62805

### Immunogen

Synthetic Peptide

### Storage

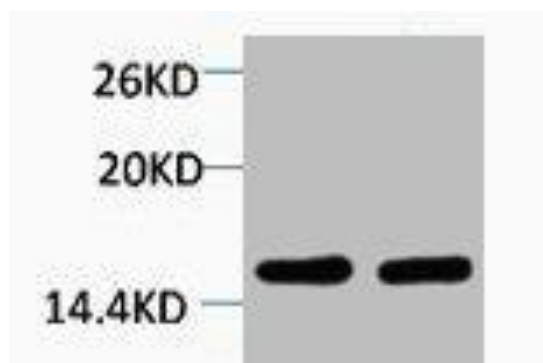
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

HIST1H4A; H4/A; H4FA; HIST1H4B; H4/I; H4FI; HIST1H4C; H4/G; H4FG; HIST1H4D; H4/B; H4FB; HIST1H4E; H4/J; H4FJ; HIST1H4F; H4/C; H4FC; HIST1H4H; H4/H; H4FH; HIST1H4I; H4/M; H4FM; HIST1H4J; H4/E; H4FE; HIST1H4K; H4/D; H4FD; HIST1H4L; H4/K; H4FK; HIST2H4A; H4/

### Application

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



Western blot analysis of 1) HeLa and 2) 3T3 at a 1:2000 dilution using the Histone H4K59me2 (H4K59 Dimethyl) Polyclonal Antibody.