

Histone H3K27me2 (H3K27 Dimethyl) Monoclonal Antibody [3B12]

(Catalog # A68392)

Description

Histone H3K27me2 (H3K27 Dimethyl) Monoclonal Antibody [3B12]. Unconjugated. Raised in: Mouse.

Formulation

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

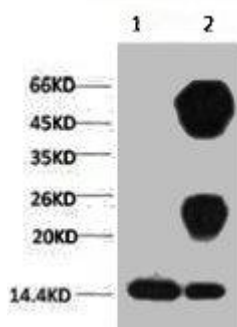
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

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/

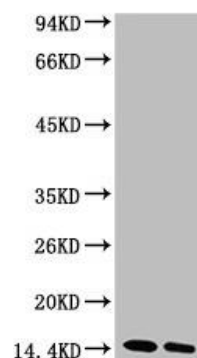
Application

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



Western blot

All lanes: FAM63A Polyclonal Antibody at 8ug/ml
Lane 1: Mouse liver tissue



Western blot analysis of HeLa, diluted at 1) 1:1000 2) 1:3000