

Histone H1K25me2 (H1K25 Dimethyl) Polyclonal Antibody

(Catalog #A68343)

Description

Histone H1K25me2 (H1K25 Dimethyl) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q02539/P16403/P16402/P10412

Purification

Affinity Purified

Immunogen

Synthetic Peptide

Storage

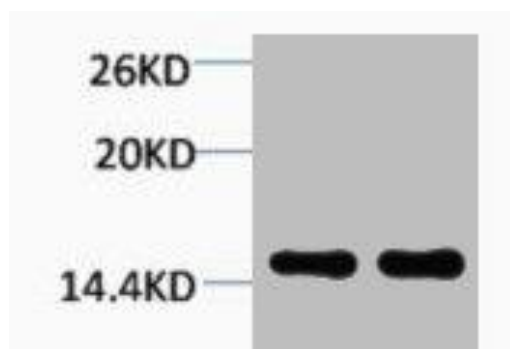
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

HIST1H1A; H1F1; Histone H1.1; Histone H1a; HIST1H1C; H1F2; Histone H1.2; Histone H1c; Histone H1d; Histone H1s-1; HIST1H1D; H1F3; Histone H1.3; Histone H1c; Histone H1s-2; HIST1H1E; H1F4; Histone H1.4; Histone H1b; Histone H1s-4

Application

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



Western blot analysis of 1) HeLa, 2) 3T3, diluted at 1:2000.