
Histone H2BK43me2 (H2BK43 Dimethyl) Polyclonal Antibody

(Catalog #A68346)

Description

Histone H2BK43me2 (H2BK43 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

Q96A08/P33778/P62807

Purification

Affinity Purified

Immunogen

Synthetic Peptide

Storage

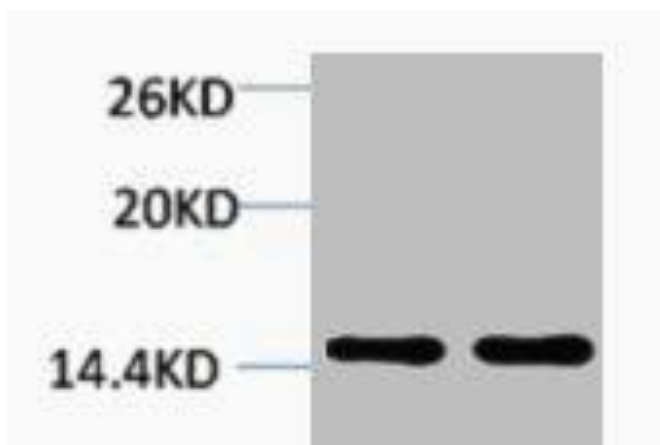
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

HIST1H2BA; TSH2B; Histone H2B type 1-A; Histone H2B, testis; Testis-specific histone H2B; HIST1H2BB; H2BFF; Histone H2B type 1-B; Histone H2B.1; Histone H2B.f; H2B/f; HIST1H2BC; H2BFL; HIST1H2BE; H2BFH; HIST1H2BF; H2BFG; HIST1H2BG; H2BFA; HIST1H2BI; H2BFK

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.