
Histone H2BK126ac (Acetyl H2BK126) Polyclonal Antibody

(Catalog #A68380)

Description

Histone H2BK126ac (Acetyl H2BK126) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

PBS with 50% glycerol, 0.5% BSA* and 0.02% sodium azide.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q96A08

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.

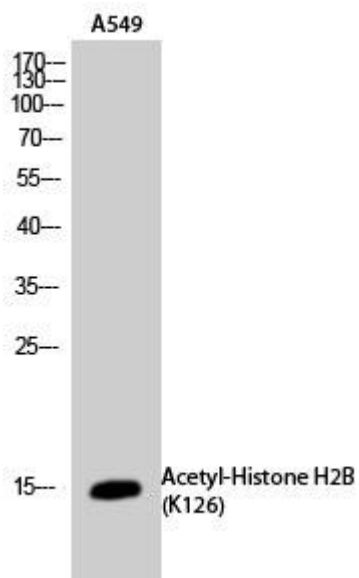
Alternative Names

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

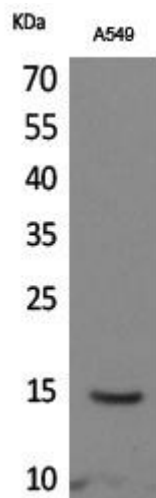
Application

WB, ELISA; Recommended dilution: WB:1:500-1:2000, ELISA:1:20000

**The BSA is derived from animal or animal-derived material of negligible amounts. The animal-derived material is subject to heat treatment at a temperature higher than 65°C for at least three hours and acid treatment with pH value of less than 5 for at least three hours, thereby being free of Bovine Spongiform Encephalopathy (BSE) and Transmissible Spongiform Encephalopathy (TSE). Additionally, as all BSA-containing products are strictly intended for research purposes and not for diagnostic or therapeutic use, they are not subject to certification authority oversight.*



Western blot analysis of A549 cells using the Histone H2BK126ac (Acetyl H2BK126) Polyclonal Antibody.



Western blot analysis of A549 cells using the Histone H2BK126ac (Acetyl H2BK126) Polyclonal Antibody.