

Histone H2B Polyclonal Antibody

(Catalog #A68383)

Description

Histone H2B Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P62807/P06899/Q16778/P57053

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.

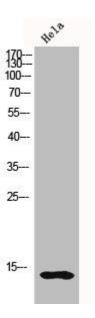
Alternative Names

HIST1H2BC; H2BFL; HIST1H2BE; H2BFH; HIST1H2BF; H2BFG; HIST1H2BG; H2BFA; HIST1H2BI; H2BFK; Histone H2B type 1-C/E/F/G/I; Histone H2B.1 A; Histone H2B.a; H2B/a; Histone H2B.g; H2B/g; Histone H2B.h; H2B/h; Histone H2B.k; H2B/k; Histone H2B.I

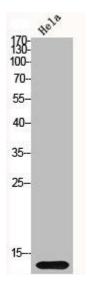
Application

WB, IHC, IF, ELISA; Recommended dilution: WB:1:500-1:2000, IHC:1:100-1:300, IF:1:200-1:1000, 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 HELA cells using Histone H2B Polyclonal Antibody



Western Blot analysis of HELA cells using Histone H2B Polyclonal Antibody