

Histone H2B antibody (mAb)

Catalog No: 61037

Clone: 5HH2-2A8

Isotype: IgG1

Application(s): WB

Reactivity: Human

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Quantity: 100 µg

Molecular Weight: 15 kDa

Background: Histone H2B is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 147 base pairs of DNA wrapped around an octamer of core histone proteins (two each of Histone H2A, Histone H2B, Histone H3 and Histone H4). Histone H1 is a linker histone, present at the interface between the nucleosome core and DNA entry/exit points. Histone H1 is responsible for establishing higher-order chromatin structure. Chromatin is subject to a variety of chemical modifications, including post-translational modifications of the histone proteins and the methylation of cytosine residues in the DNA. Reported histone modifications include acetylation, methylation, phosphorylation, ubiquitylation, glycosylation, ADP-ribosylation, carbonylation and SUMOylation; these modifications play a major role in regulating gene expression.

Immunogen: This Histone H2B antibody was raised against a peptide from the C-terminal region of human Histone H2B.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Validated Applications:

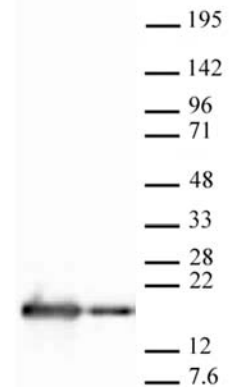
WB: 0.1 - 1 µg/ml dilution

Published Applications:

WB

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Histone H2B mAb tested by Western blot.

Detection of Histone H2B by Western blot.

Lane 1: Nuclear extract of HeLa cells (20 µg).

Lane 2: 500 ng of recombinant human

Histone H2B. Both probed with the Histone

H2B antibody at 0.1 µg/ml.