

Histone H2BK120ub1 antibody (mAb)

Catalog Nos: 39623, 39624

RRID: AB_2793279

Clone: 56

Application(s): ChIP, WB

Reactivity: Human, Wide Range Predicted

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Isotype: IgG2a

Concentration: 1 µg/µl

Molecular Weight: 26 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; it 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; they play a major role in regulating gene expression.

Histone H2B is ubiquitylated at lysine 120 in vertebrates. H2B ubiquityl Lys120 is found preferentially associated with the transcribed regions of highly expressed genes. It is possible that H2B ubiquitylation is associated with transcriptional elongation.

Immunogen: This Histone H2B ubiquityl Lys120 antibody was raised against a branched peptide corresponding the the ubiquitin conjugation site at Lys120 of human Histone H2B.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

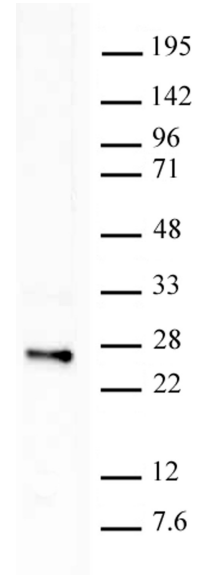
Applications Validated by Active Motif:

WB*: 0.5 - 2 µg/ml dilution

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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



Histone H2B ubiquityl Lys120 antibody tested by Western blot.

HeLa nuclear extract (20 µg per lane) probed with Histone H2B ubiquityl Lys120 antibody (1 µg/ml).