

Histone H2AK5ac antibody (pAb)

Catalog Nos: 39107, 39108

RRID: AB_2793157

Isotype: Serum

Application(s): ChIP, DB, IF, WB

Reactivity: Human, Wide Range Predicted

Volumes: 200 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 14 kDa

Background: Histone H2A 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.

Lysine N-ε-acetylation is a dynamic, reversible and tightly regulated protein and histone modification that plays a major role in chromatin remodeling and in the regulation of gene expression in various cellular functions. Histone H2A Lys5 is preferentially acetylated by TIP60 and deacetylated by HDAC3.

Immunogen: This Histone H2A acetyl Lys5 antibody was raised against a peptide including acetyl-lysine 5 of human histone H2A.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

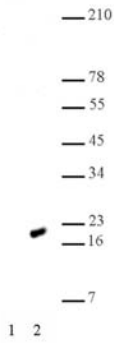
WB*: 1:500 - 1:2,000 dilution

IF: 1:500 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.

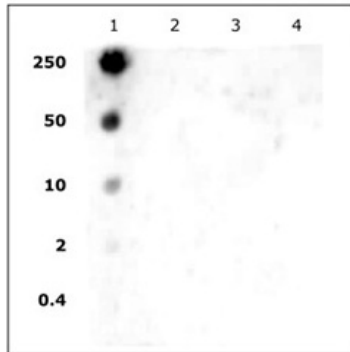


Histone H2A acetyl Lys5 pAb tested by Western blot

HeLa acid extract (5 µg per lane) was probed with Histone H2A acetyl Lys5 polyclonal antibody at a 1:2,000 dilution.

Lane 1: No treatment.

Lane 2: Cells treated with sodium butyrate.



Specificity data:

Dot blot analysis was used to confirm the specificity of Histone H2A acetyl Lys5 pAb for acetyl Lys5 of Histone H2A. Modified and unmodified peptides were spotted onto PVDF and probed with the antibody at a 1:2,000 dilution. The amount of peptide spotted (in picomoles) is indicated next to each row.

Lane 1: Peptide acetylated at Lys5 of H2A.

Lane 2: Unmodified Lys5 peptide.

Lane 3: Peptide acetylated at Lys9 of H2A.

Lane 4: Unmodified Lys9 peptide.

Detection of Histone H2AK5ac by immunofluorescence.

U2OS cells were stained with H2AK5ac antibody at a dilution of 1:500. Left panel: DAPI. Middle panel: H2AK5ac antibody staining. Right panel: merge.

