

Histone H3K18ac antibody (pAb)

Catalog Nos: 39755, 39756

RRID: AB_2714186

Isotype: IgG

Application(s): ChIP, ChIP-Seq, DB, ICC, IF, WB

Reactivity: Human, Wide Range Predicted

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Rabbit

Concentration: 1 µg/µl

Molecular Weight: 17 kDa

Background: Histone H3 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). 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. Acetylation of histones is linked to a number of specific processes including transcriptional regulation and genomic organization.

Immunogen: This Histone H3 acetyl Lys18 antibody was raised against a peptide including acetyl-lysine 18 of histone H3.

Buffer: Purified IgG in PBS (pH 7.5) with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic. For your convenience, an unpurified serum version (Catalog No. 39587) of this antibody is also available.

Application Notes:

Applications Validated by Active Motif:

ChIP: 10 µg per ChIP

ChIP-Seq: 4 µg each

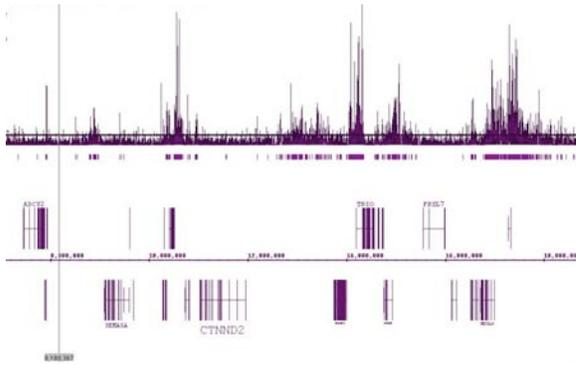
ICC/IF: 1 µg/ml dilution

WB: 0.5 - 2 µg/ml dilution

DB: 0.1 - 0.5 µg/ml dilution

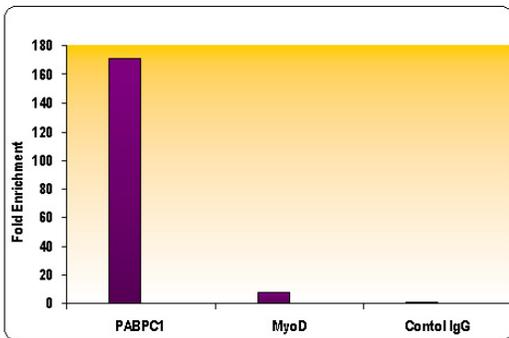
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



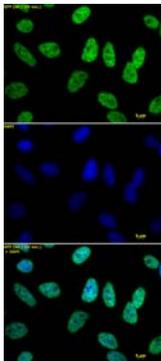
Histone H3 acetyl Lys18 antibody tested by ChIP-chip.

ChIP was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with chromatin from 2 million HeLa cells. ChIP DNA was amplified by WGA, labeled and hybridized to a human tiling array. The image shows a 12 million bp view on chromosome 5 with long stretches of H3K18Ac enrichment



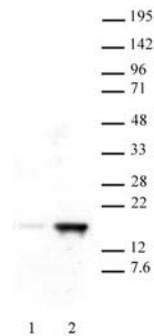
Histone H3 acetyl Lys18 antibody tested by ChIP.

Chromatin IP performed using the ChIP-IT[®] Express Kit (Catalog No. 53008) and HeLa Chromatin (1.5 x 10⁶ cell equivalents per ChIP) using 10 µg of Histone H3 acetyl Lys18 antibody or the equivalent amount of rabbit IgG as a negative control. Real time, quantitative PCR (RT-qPCR) was performed on DNA purified from each of the ChIP reactions using a primer pair specific for the indicated gene. Data are presented as Fold Enrichment of the ChIP antibody signal versus the negative control IgG using the ddCT method.



Histone H3 acetyl Lys18 antibody tested by immunofluorescence.

Detection of Histone H3 acetyl Lys18 by immunofluorescence. HeLa cells were stained with Histone H3 acetyl Lys18 antibody at a dilution of 1 µg/ml. Top panel: Histone H3 acetyl Lys18 antibody staining. Middle panel: DAPI. Bottom panel: merge.



Histone H3 acetyl Lys18 antibody tested by Western blot.

HeLa cell extract (20 µg per lane) probed with Histone H3 acetyl Lys18 antibody (0.5 µg/ml).
Lane 1: Untreated cells.
Lane 2: Cells treated with Sodium Butyrate..

Histone H3 acetyl Lys18 antibody tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of Histone H3 acetyl Lys18 antibody for acetyl Lys18 histone H3. Acetylated peptides corresponding to the immunogen and related peptides were spotted onto PVDF and probed with the antibody at a dilution of 0.1 µg/ml. The amount of peptide (picomoles) spotted is indicated next to each row.

Lane 1: acetyl-Lys4 peptide. Lane 2: unmodified Lys4 peptide. Lane 3: acetyl-Lys18 peptide. Lane 4: unmodified Lys18 peptide. Lane 5: acetyl-Lys9 peptide. Lane 6: acetyl-Lys14 peptide. Lane 7: acetyl-Lys18 peptide 2. Lane 8: acetyl-Lys23 peptide. Lane 9: acetyl-Lys27 peptide.

