

Histone H3K9me0 antibody (mAb)

Catalog Nos: 61399, 61400

RRID: AB_2793620

Clone: 9B1-2G6

Isotype: IgG1

Application(s): WB

Reactivity: Human, Wide Range Predicted

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 17 kDa

Background: Histone H3 is one of the core components of the nucleosome, the basic building block of chromatin. Histones are 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.

Methylation of lysines occur as four states: unmethylated (me0), monomethyl (me1), dimethyl (me2) and trimethyl (me3). The me0 state of lysine is recognized as biologically relevant and a number of proteins containing PhD fingers, ADD and WD40 domains are known to associate with unmodified lysines.

Immunogen: This antibody was raised against a synthetic branched peptide corresponding to amino acids surrounding Lys9 of human Histone H3.

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

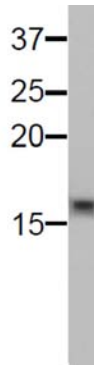
Application Notes:

Applications Validated by Active Motif:

WB: 0.5 - 2 µg/ml dilution

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.

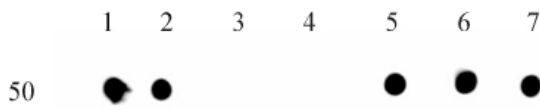


Histone H3K9me0 mAb tested by Western blot.

Histone H3K9me0 detection by Western blot. The analysis was performed using HeLa acid extract.

Histone H3K9me0 mAb tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of Histone H3K9me0 mAb for detection of unmethylated Lys9 of Histone H3. 50 pmol of recombinant Histone H3 proteins containing various modifications were spotted onto PVDF and probed with the antibody at 1 µg/mL.



Lane 1: Histone H3 unmodified.

Lane 2: Histone H3K4me3.

Lane 3: Histone H3K9me3.

Lane 4: Histone H3 K9ac.

Lane 5: Histone H3 K14ac.

Lane 6: Histone H3K4ac.

Lane 7: Histone H3 S10ph.