

Histone H3K79me2 antibody (pAb)

Catalog Nos: 39923, 39924

RRID: AB_2793395

Isotype: IgG

Application(s): ChIP, DB, 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.

Lysine 79 of histone H3 can be mono-, di- or trimethylated by Dot1 methylase; methylation at this residue acts as a marker of inactive chromatin regions that is critical for transcriptional silencing, and it is thought that silencing proteins such as Sir3 function by blocking Dot1 methylation.

Immunogen: This Histone H3 dimethyl Lys79 antibody was raised against a peptide including dimethyl-lysine 79 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. 39143) of this antibody is also available.

Application Notes:

Applications Validated by Active Motif:

ChIP: 10 µg per ChIP

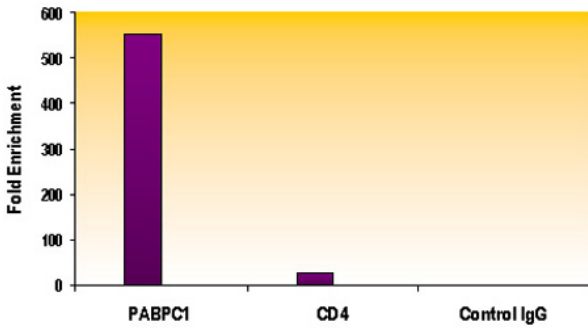
WB*: 1 - 2 µg/ml dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

*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.

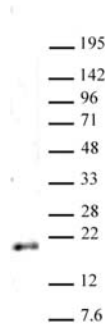
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product 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 H3 dimethyl Lys79 antibody tested by ChIP analysis.

Chromatin IP performed using the ChIP-IT[®] Express Kit (Catalog No. 53008) and HeLa Chromatin (1.5×10^6 cell equivalents per ChIP) using 10 μ g of Histone H3 dimethyl Lys79 pAb 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 dimethyl Lys79 antibody (pAb) tested by Western blot.

Nuclear extract of HeLa cells (20 μ g) probed with Histone H3 dimethyl Lys79 antibody (1 μ g/ml dilution).



Histone H3 dimethyl Lys79 antibody (pAb) tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of Histone H3 dimethyl Lys79 antibody for dimethyl-Lys79 of histone H3. Peptides corresponding to the region around lysine 79 of histone H3 were spotted onto PVDF and probed with Histone H3 dimethyl Lys79 antibody at a dilution of 0.5 μ g/ml. The amount of peptide (in picomoles) spotted is indicated next to each row.

Top panel: Lane 1: unmodified Lys4. Lane 2: monomethyl Lys4. Lane 3: dimethyl Lys4. Lane 4: trimethyl Lys4. Lane 5: unmodified Lys9. Lane 6: monomethyl Lys9. Lane 7: dimethyl Lys9. Lane 8: trimethyl Lys9. Lane 9: unmodified Lys79. Lane 10: monomethyl Lys79. Lane 11: dimethyl Lys79. Lane 12: trimethyl Lys79.

Bottom panel: Lane 1: Unmodified Lys23. Lane 2: Monomethyl Lys23. Lane 3: Dimethyl Lys23. Lane 4: Trimethyl Lys23. Lane 5: unmodified Lys27. Lane 6: monomethyl Lys27. Lane 7: dimethyl Lys27. Lane 8: trimethyl Lys27. Lane 9: unmodified Lys36. Lane 10: monomethyl