

Histone H3K23me2 antibody (pAb)

Catalog Nos: 39653, 39654

RRID: AB_2793291

Isotype: Serum

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

Reactivity: Human, Wide Range Predicted

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

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). Histone H1 is a linker histone, present at the interface between the nucleosome core and DNA entry/exit points. Histone H1 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; these modifications play a major role in regulating gene expression. The methylation of histones can occur on two different residues: arginine or lysine. Histone methylation can be associated with transcriptional activation or repression, depending on the methylated residue.

Immunogen: This Histone H3 dimethyl Lys23 antibody was raised against a peptide containing dimethyl-Lys23 of human histone H3.

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

Application Notes:

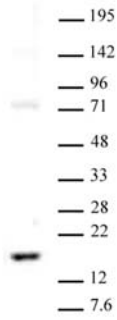
Applications Validated by Active Motif:

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

DB: 1:500 - 1:2,000 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 dimethyl Lys23 pAb tested by Western blot.

HEK293 nuclear extract (25 µg) was probed with Histone H3 dimethyl Lys23 pAb (1:500 dilution).

Histone H3 dimethyl Lys23 pAb tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of Histone H3 dimethyl Lys23 pAb for dimethyl-Lys23 of histone H3. Peptides corresponding to regions around major sites of histone H3 methylation were spotted onto PVDF and probed with Histone H3 dimethyl Lys23 pAb at a dilution of 1: 2,000. The amount of peptide (in picomoles) spotted is indicated next to each row.

Top Panel: Lane 1: dimethyl Lys14. Lane 2: monomethyl Lys18. Lane 3: dimethyl Lys18. Lane 4: trimethyl Lys18. Lane 5: unmodified Lys23. Lane 6: monomethyl Lys23. Lane 7: dimethyl Lys23. Lane 8: trimethyl Lys23. Lane 9: unmodified Lys27. Lane 10: monomethyl Lys27. Lane 11: dimethyl Lys27. Lane 12: trimethyl Lys27. Bottom 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.

