

Histone H3R17me2a (asymmetric) antibody (pAb)

Catalog Nos: 39709, 39710

RRID: AB_2793315

Isotype: IgG

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

Reactivity: Human, Wide Range Predicted

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

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. The methylation of histones can occur on two different residues: arginine or lysine.

Immunogen: This Histone H3 dimethyl Arg17 asymmetric antibody was raised against a peptide containing dimethyl-Arg17 (asymmetric) of human histone H3.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

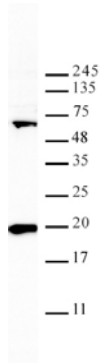
Applications Validated by Active Motif:

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

*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.

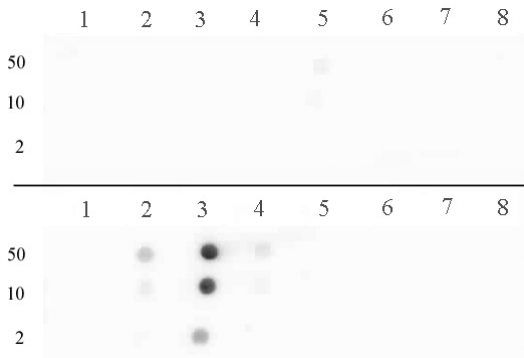
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product 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 Arg17 asymmetric antibody (pAb) tested by Western blot.

HeLa nuclear extract (20 µg) probed with Histone H3 dimethyl Arg17 asymmetric antibody at a 1:500 dilution.



Histone H3 dimethyl Arg17 asymmetric pAb tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of Histone H3 dimethyl Arg17 asymmetric pAb for dimethyl Arg17 of histone H3. Peptides corresponding to the immunogen and related peptides were spotted onto PVDF and probed with Histone H3 dimethyl Arg17 asymmetric pAb at a dilution of 1:30,000. The amount of peptide (picomoles) spotted is indicated next to each row.

Top panel Lane 1: unmodified Arg2. Lane 2: monomethyl-Arg2 H3. Lane 3: dimethyl-Arg2 H3 (sym). Lane 4: dimethyl-Arg2 H3 (asym). Lane 5: unmodified Arg8. Lane 6: monomethyl-Arg8 H3. Lane 7: dimethyl-Arg8 H3 (sym). Lane 8: dimethyl-Arg8 H3 (asym).

Bottom panel Lane 1: unmodified Arg17. Lane 2: monomethyl-Arg17 H3. Lane 3: dimethyl-Arg17 H3 (asym). Lane 4: dimethyl-Arg17 H3 (sym). Lane 5: unmodified Arg26. Lane 6: monomethyl-Arg26 H3. Lane 7: dimethyl-Arg26 H3 (asym). Lane 8: dimethyl-Arg26 H3 (sym).