

Histone H3R2me2s (symmetric) antibody (pAb)

Catalog No: 39703

RRID: AB_2793312

Isotype: IgG

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

Reactivity: Human, Wide Range Predicted

Volume: 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 Arg2 (H3R2me2s) antibody was raised against a peptide containing dimethyl-Arg2 (symmetric) of human histone H3.

Buffer: Purified IgG in PBS with 50% glycerol and 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

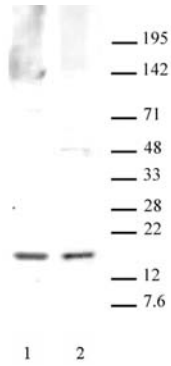
WB*: 1:500 dilution

DB: 1:10,00 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.

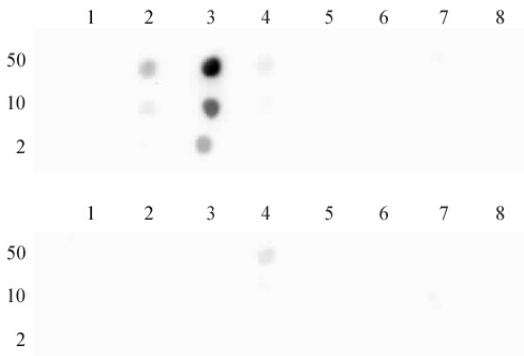
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product 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 H3R2me2s (symmetric) antibody (pAb) tested by Western blot.

Western blot performed using Histone H3R2me2s (symmetric) antibody (pAb) at a dilution of 1:500. Lane 1: Nuclear extract from HEK293 cells (25 µg). Lane 2: Nuclear extract from HeLa cells (20 µg).



Histone H3R2me2s (symmetric) antibody (pAb) tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of Histone H3R2me2s (symmetric) antibody (pAb) for dimethyl-arginine 2 of histone H3. Peptides corresponding to the immunogen and related peptides were spotted onto PVDF and probed with Histone H3 dimethyl Arg2 symmetric pAb at 1:10,000. 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 (symmetric). Lane 4: dimethyl-Arg2 H3 (asymmetric). Lane 5: unmodified Arg8. Lane 6: monomethyl-Arg8 H3. Lane 7: dimethyl-Arg8 H3 (symmetric). Lane 8: dimethyl-Arg8 H3 (asymmetric). Bottom Panel: Lane 1: unmodified Arg17 H3. Lane 2: monomethyl-Arg17. Lane 3: dimethyl-Arg17 H3 (symmetric). Lane 4: dimethyl-Arg17 H3 (asymmetric). Lane 5: unmodified Arg26. Lane 6: monomethyl-Arg26 H3. Lane 7: dimethyl-Arg26 H3 (symmetric). Lane 8: dimethyl-Arg26 H3 (asymmetric).