

Histone H3R8me1 antibody (pAb)

Catalog Nos: 39673, 39674

RRID: AB_2793299

Isotype: Serum

Application(s): DB, 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 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. Histone H3 Arg8, methylated by PRMT enzymes, is involved in nuclear-receptor-mediated transcriptional activation.

Immunogen: This Histone H3 monomethyl Arg8 antibody was raised against a peptide containing monomethyl-Arg8 of human histone H3.

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

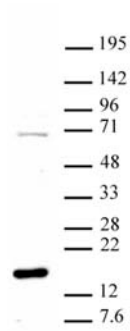
Application Notes:

Applications Validated by Active Motif:

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

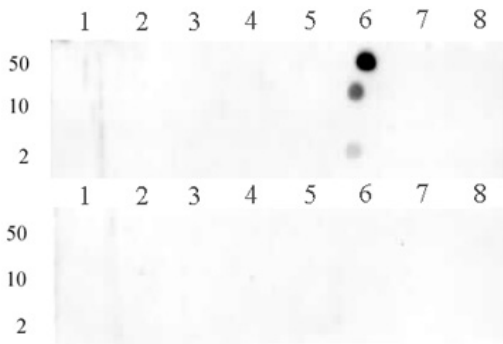
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product 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 monomethyl Arg8 pAb tested by Western blot.

HeLa cell nuclear extract (20 µg) probed with Histone H3 monomethyl Arg8 pAb (1:500 dilution).



Histone H3 monomethyl Arg8 pAb tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of Histone H3 monomethyl Arg8 pAb for monomethyl-arginine 8 of histone H3. Peptides corresponding to the immunogen and related peptides were spotted onto PVDF and probed with Histone H3 monomethyl Arg8 pAb at 1:1,000. The amount of peptide (picomoles) spotted is indicated next to each row.

Top Panel:

- Lane 1: unmodified Arg2 peptide.
- Lane 2: monomethyl-Arg2 H3 peptide.
- Lane 3: dimethyl-Arg2 H3 (symmetric) peptide.
- Lane 4: dimethyl-Arg2 H3 (asymmetric) peptide.
- Lane 5: unmodified Arg8 peptide.