

Histone H3K122me1 antibody (pAb)

Catalog Nos: 39367, 39368

RRID: AB_2793230

Isotype: IgG

Application(s): DB, WB

Reactivity: Human, Wide Range Predicted

Volumes: 200 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 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.

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 monomethyl Lys122 antibody was raised against a peptide containing monomethyl-lysine 122 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.

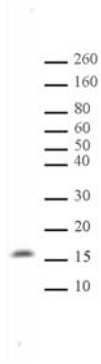
Application Notes:

Applications Validated by Active Motif:

WB: 1:250 - 1:1,000 dilution

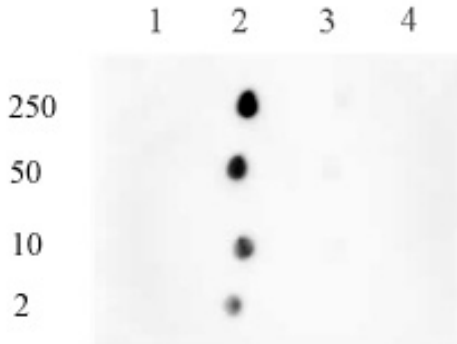
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product 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 Lys122 pAb tested by Western blot.

Nuclear extract of HeLa cells (20 µg) was probed with Histone H3 monomethyl Lys122 pAb (1:500 dilution).



Histone H3 monomethyl Lys122 pAb tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of 39367 for monomethyl Lys122 histone H3. Methylated peptides corresponding to the immunogen were spotted onto PVDF and probed with 39367 at 1:5,000. The amount of peptide (picomoles) spotted is indicated next to each row.

- Lane 1: unmodified Lys122 H3 peptide.
- Lane 2: monomethyl-Lys122 H3 peptide.
- Lane 3: dimethyl-Lys122 H3 peptide.
- Lane 4: trimethyl-Lys122 H3 peptide.