

Histone H4K8K12K16 (pan-biotinylated) antibody (pAb)

Catalog Nos: 61393, 61394

RRID: AB_2793617

Isotype: IgG

Application(s): DB

Reactivity: Human, Wide Range Predicted

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 8 kDa

Background: Histone H4 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; it 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; they play a major role in regulating gene expression.

The biotinylation of Histone H4 on Lys8, Lys12 and Lys16 has been reported to be enriched in repeat regions and may participate in repression of transcriptionally competent genes.

Immunogen: This antibody was raised against a synthetic peptide containing biotin-Lys 12 of human Histone H4.

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

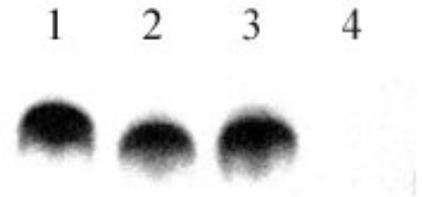
Application Notes:

Applications Validated by Active Motif:

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 H4K8K12K16 (pan-biotinylated) pAb tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of Histone H4K8K12K16 (pan-biotinylated) antibody. Modified and unmodified peptide were spotted onto the membrane and probed with the antibody.

Lane 1: Peptide biotinylated at lysine 8 of Histone H4.

Lane 2: Peptide biotinylated at lysine 12 of Histone H4.

Lane 3: Peptide biotinylated at lysine 16 of Histone H4.