

Histone H3, C-terminal antibody (pAb)

Catalog Nos: 39451, 39052, 39452

RRID: AB_2793242

Isotype: IgG

Application(s): ELISA, WB

Reactivity: Human

Volumes: 100 μ l, 50 μ l, 10 μ l

Purification: None

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). Histone H1 is a linker histone, present at the interface between the nucleosome core and DNA entry/exit points. Histone H1 is responsible for establishing higher-order chromatin structure.

Each histone contains two domains. First, a main globular domain (C-terminal) forming the core of the nucleosome is involved in histone-histone interactions and in binding to the DNA. Secondly, an N-terminal tail is subject to post-translational modifications. 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.

Immunogen: This Histone H3, C-terminal antibody was raised against a C-terminal peptide of histone H3.

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

Application Notes:

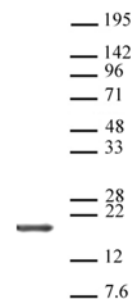
Applications Validated by Active Motif:

WB*: 1:5,000 - 1:25,000 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.

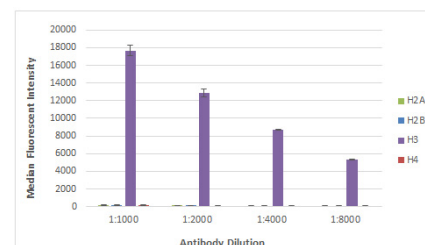
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

Analysis was performed using HeLa cell nuclear extract and Histone H3, C-terminal polyclonal antibody at 1:25,000 dilution. M.W. of the detected protein is 17 kDa.



Specificity of Histone H3, C-terminal (pAb)

Human Histone proteins were conjugated to Luminex MagPlex beads and incubated with various amounts of Histone H3, C-terminal antibody. Protein-bound antibody was detected with Anti-Rabbit IgG-PE and read in a MagPix Luminex Instrument.