

Histone H4 antibody (pAb)

Catalog Nos: 39269, 39070, 39270

RRID: AB_2636967

Isotype: Serum

Application(s): ChIP, WB

Reactivity: Human, Wide Range Predicted

Volumes: 100 µl, 50 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.

Immunogen: This Histone H4 antibody was raised against full-length recombinant *Xenopus* histone H4.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic. For your convenience, an IgG version (Catalog No. 61199) of this antibody that was purified by Protein A Chromatography is also available.

Application Notes:

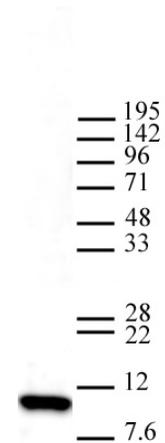
Applications Validated by Active Motif:

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

For Histone H4, we also offer AbFlex® Histone H4 Recombinant Antibody (rAb). For details, see Catalog No. 91295.

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 H4 antibody (pAb) tested by Western blot.

Detection of Histone H4 by Western blot. The analysis was performed using HeLa acid extract (20 µg) and probed with Histone H4 pAb at a 1:1,000 dilution.