

## Histone H4 antibody (pAb)

**Catalog Nos:** 61299, 61300

**RRID:** AB\_2650524

**Isotype:** IgG

**Application(s):** ChIP, ChIP-Seq, WB

**Reactivity:** Human, Wide Range Predicted

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Concentration:** 1  $\mu$ g/ $\mu$ l

**Molecular Weight:** 8 kDa

**Background:** Histone H4 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.

**Immunogen:** This Histone H4 antibody was raised against a peptide corresponding to amino acids 43- 54 of human histone H4.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

ChIP: 2 - 10  $\mu$ l per ChIP

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

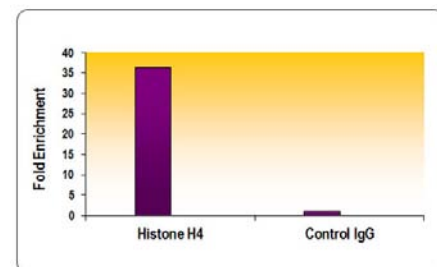
\*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western Blot.

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

**References:** Histone H4 antibody has been cited in one or more publications. To see an up-to-date listing of papers that describe its use, please go to [www.activemotif.com/catalog/details/61299](http://www.activemotif.com/catalog/details/61299).

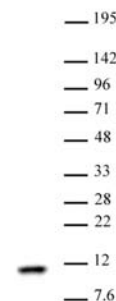
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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

Performed using HeLa Chromatin ( $1.5 \times 10^6$  cell equivalents per ChIP) and 10  $\mu$ l of Histone H4 antibody or 10  $\mu$ l of rabbit IgG as a negative control. Real time, quantitative PCR was performed on DNA purified from each of the ChIP reactions using a primer pair specific for the PABPC1 (+ 0.5) gene. Data are presented as Fold Enrichment of the ChIP antibody signal versus the negative control IgG (ddCT method).



### Histone H4 antibody (pAb) tested by Western blot

Nuclear extract of HeLa cells (30  $\mu$ g) probed with Histone H4 antibody (1:1,000).