

## Histone H3 antibody (pAb)

**Catalog Nos:** 61799, 61800

**RRID:** AB\_2793771

**Isotype:** IgG

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

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Rabbit

**Concentration:** 1 µg/µl

**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.

**Immunogen:** This antibody was raised against Recombinant Histone H3 from *Xenopus laevis*, which shares 99% identity to human H3 (134/136 aa are identical).

**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-Seq: 4 µg per ChIP

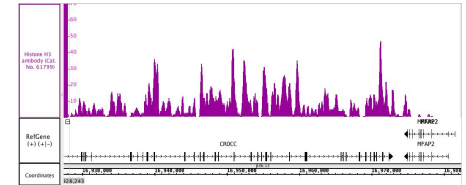
WB\*: 0.1 – 1 µg/ml

\*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 H3, we also offer AbFlex® Histone H3 Recombinant Antibody (rAb). For details, see Catalog No. 91297.

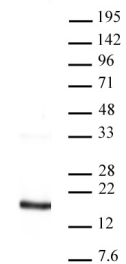
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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

Chromatin immunoprecipitation (ChIP) was performed using the [ChIP-IT® High Sensitivity Kit](#) (Cat. No. 53040) with 30 µg of chromatin from Farage human lymph-node derived B cells and 4 µl Histone H3 antibody. ChIP DNA was sequenced on the Illumina HiSeq and 20 million sequence tags were mapped to identify Histone H3 binding sites.



### Histone H3 antibody tested by Western blot

20 µg HeLa cell nuclear extract was run on an SDS-PAGE gel and probed with Histone H3 antibody (pAb) at 0.5 µg/ml.

MW: 17 kDa