

## Histone H3K14ac antibody (mAb)

---

**Catalog No:** 61433

**RRID:** AB\_2793634

**Clone:** 13HH3-1A5

**Isotype:** IgG1, k

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

**Reactivity:** Human, Mouse, Wide Range Predicted

**Quantity:** 100 µg

**Purification:** Protein A Chromatography

**Host:** Mouse

**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. Acetylation of histones is linked to a number of specific processes including transcriptional regulation and genomic organization.

**Immunogen:** This antibody was raised against a peptide corresponding to amino acids 10-18 including acetyl-lysine 14 of human Histone H3.

**Buffer:** Purified IgG in 60 mM Sodium Citrate, 105 mM Tris-HCl, 30% glycerol, and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

**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.