

# PRODUCT INFORMATION



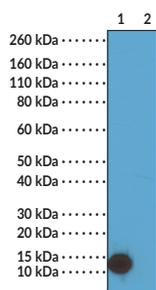
## Histone H4K5Ac Monoclonal Antibody (RM199)

Item No. 32156

### Overview and Properties

**Contents:** This vial contains 100 µg of protein A-affinity purified monoclonal antibody.  
**Synonym:** Acetylated Histone H4 Lysine 5  
**Immunogen:** Peptide corresponding to H4K5Ac  
**Cross Reactivity:** (+) H4K5Ac; (-) Unmodified H4K5, H4K8Ac, H4K12Ac, H4K16Ac, H4K20Ac, H4K31Ac, H4K91Ac  
**Species Reactivity:** (+) Vertebrates  
**Form:** Liquid  
**Storage:** -20°C (as supplied)  
**Stability:** ≥1 year  
**Storage Buffer:** PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide  
**Concentration:** 1 mg/ml  
**Clone:** RM199  
**Host:** Rabbit  
**Isotype:** IgG  
**Applications:** ELISA, immunocytochemistry (ICC), multiplex-based assays, and Western blot (WB); the recommended starting concentration is 0.5-2 and 0.05-0.2 µg/ml for ICC and multiplex-based assays, respectively, and 0.2-1 µg/ml for ELISA and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

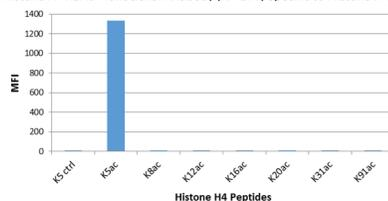
### Images



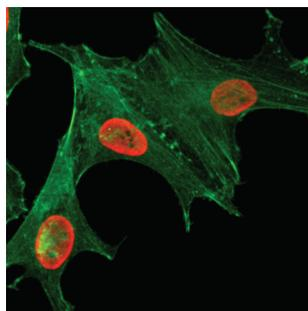
Lane 1: Acid extracts of HeLa cells  
Lane 2: Recombinant histone H4

WB of acid extracts of HeLa cells and recombinant histone H4. Using Histone H4K5Ac Monoclonal Antibody (RM199) at 0.2 µg/ml showed a band of H4K5Ac in HeLa cells.

Histone H4K5Ac Monoclonal Antibody (RM199) Specific to Histone H4K5Ac



Histone H4K5Ac monoclonal antibody (RM199) specifically reacts to H4K5Ac. No cross reactivity with unmodified H4K5, H4K8Ac, H4K12Ac, H4K16Ac, H4K20Ac, H4K31Ac, or H4K91Ac.



Immunocytochemistry of HeLa Cells Treated with Sodium Butyrate using Histone H4K5Ac Monoclonal Antibody (RM199) (red). Actin filaments have been labeled with fluorescein phalloidin (green).

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/26/2024

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

Histone H4 is one of four core histone proteins that are involved in the organization of DNA into chromatin.<sup>1</sup> Histones are globular proteins with unstructured N-terminal tails and are subject to a variety of posttranslational modifications, such as methylation, acetylation, phosphorylation, and citrullination, that can influence chromatin structure and regulate gene transcription.<sup>1,2</sup> Acetylation of histone H4 at lysine 5 (H4K5Ac) is enriched in promoter regions and coding sequences in mouse hippocampus and is associated with active chromatin.<sup>3,4</sup> Hippocampal levels of H4K5Ac increase following contextual fear conditioning in mice.<sup>3</sup> Levels of H4K5Ac are elevated in postmortem brain tissue isolated from fetuses with spina bifida compared with healthy controls.<sup>5</sup> Global levels of H4K5Ac are decreased in blast cells from patients with acute myeloid leukemia compared with healthy individuals, and low levels of H4K5Ac are associated with poor prognosis in these patients.<sup>6</sup> Cayman's Histone H4K5Ac Monoclonal Antibody (RM199) can be used for ELISA, immunocytochemistry (ICC), multiplex-based assay, and Western blot (WB) applications.

## References

---

1. Wang, Y., Li, M., Stadler, S., *et al.* Histone hypercitrullination mediates chromatin decondensation and neutrophil extracellular trap formation. *J. Cell Biol.* **184**(2), 205-213 (2009).
2. Hyun, K., Jeon, J., Park, K., *et al.* Writing, erasing and reading histone lysine methylations. *Exp. Mol. Med.* **49**(4), e324 (2017).
3. Park, C.S., Rehrauer, H., and Mansuy, I.M. Genome-wide analysis of H4K5 acetylation associated with fear memory in mice. *BMC Genomics* **14**, 539 (2013).
4. Zhao, R., Nakamura, T., Fu, Y., *et al.* Gene bookmarking accelerates the kinetics of post-mitotic transcriptional re-activation. *Nat. Cell Biol.* **13**(11), 1295-304 (2011).
5. Li, D., Wan, C., Bai, B., *et al.* Identification of histone acetylation markers in human fetal brains and increased H4K5ac expression in neural tube defects. *Mol. Genet. Genomic Med.* **7**(12), e1002 (2019).
6. Sauer, T., Arteaga, M.F., Isken, F., *et al.* MYST2 acetyltransferase expression and Histone H4 Lysine acetylation are suppressed in AML. *Exp. Hematol.* **43**(9), 794-802 (2015).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM