

# PRODUCT INFORMATION



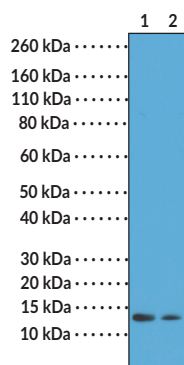
## Histone H4K12Ac Monoclonal Antibody (RM202)

Item No. 32162

### Overview and Properties

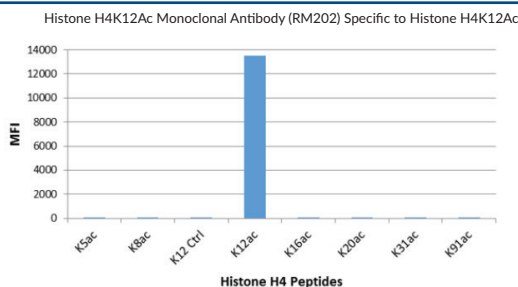
**Contents:** This vial contains 100 µg of protein A-affinity purified monoclonal antibody.  
**Synonym:** Acetylated Histone H4 Lysine 12  
**Immunogen:** Peptide corresponding to H4K12Ac  
**Cross Reactivity:** (+) H4K12Ac; (-) Unmodified H4K12, H4K5Ac, H4K8Ac, H4K16Ac, H4K20Ac, H4K31Ac, H4K91Ac  
**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:** 1.0 mg/ml  
**Clone:** RM202  
**Host:** Rabbit  
**Isotype:** IgG  
**Applications:** Chromatin immunoprecipitation (ChIP), ELISA, immunocytochemistry (ICC), multiplex-based assays, and Western blot (WB); the recommended starting concentration is 1-5 µg/ml for ChIP, 0.5-1 µg/ml for ELISA, and 0.5-2 µg/ml for ICC, multiplex-based assays, and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images

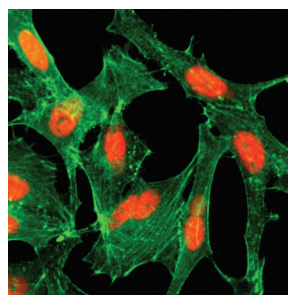


Lane 1: Acid extracts of HeLa cells treated  
Lane 2: Acid extracts of HeLa cells untreated

WB of acid extracts from HeLa cells treated with sodium butyrate or HeLa cells untreated using Histone H4K12Ac Monoclonal Antibody (RM202) at a concentration of 0.5 µg/ml.



Histone H4K12Ac Monoclonal Antibody (RM202) specifically reacts to acetylated H4K12Ac. No cross reactivity with unmodified H4K16, H4K5Ac, H4K8Ac, H4K16Ac, H4K20Ac, H4K31Ac, or H4K91Ac.



Immunocytochemical labeling of HeLa cells treated with sodium butyrate, using Histone H4K12Ac Monoclonal Antibody (RM202) (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, 02/28/2024

**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 post-translational 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 12 (H4K12Ac) is associated with chromatin relaxation and gene transcription.<sup>3</sup> Levels of H4K12Ac are elevated in two transgenic mouse models of Alzheimer's disease, as well as in monocytes isolated from patients with mild cognitive impairment and Alzheimer's disease. H4K12Ac levels are also increased in response to chronic alcohol exposure and positively correlate with the release of the pro-inflammatory cytokine chemokine (C-C motif) ligand 8 (CCL8) in monocyte-derived dendritic cells.<sup>4</sup> Cayman's H4K12Ac Monoclonal Antibody (RM202) can be used for chromatin immunoprecipitation (ChIP), 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. Plagg, B., Ehrlich, D., Kniewallner, K.M., *et al.* Increased acetylation of histone H4 at lysine 12 (H4K12) in monocytes of transgenic Alzheimer's mice and in human patients. *Curr. Alzheimer Res.* **12**(8), 752-760 (2015).
4. Parira, T., Figueroa, F., Laverde, A., *et al.* Novel detection of post-translational modifications in human monocyte-derived dendritic cells after chronic alcohol exposure: Role of inflammation regulator H4K12ac. *Sci. Rep.* **7**(1), 11236 (2017).

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