

# PRODUCT INFORMATION



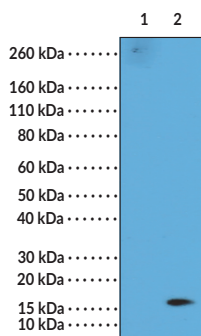
## Histone H3K36Me2 Monoclonal Antibody (RM141)

Item No. 32133

### Overview and Properties

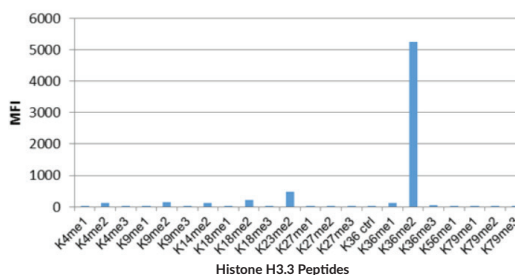
**Contents:** This vial contains 100 µg of protein A-affinity purified monoclonal antibody.  
**Synonym:** Dimethylated Histone H3 Lysine 36  
**Immunogen:** Peptide corresponding to H3K36Me2  
**Cross Reactivity:** (+) H3K36Me2; (-) Unmodified H3K36, H3K36Me1, H3K36Me3, H3K4Me1, H3K4Me2, H3K4Me3, H3K9Me1, H3K9Me2, H3K9Me3, H3K14Me2, H3K18Me1, H3K18Me2, H3K23Me1, H3K23Me2, H3K27Me1, H3K27Me2, H3K27Me3, H3K56Me1, H3K79Me1, H3K79Me2, H3K79Me3  
**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:** RM141  
**Host:** Rabbit  
**Isotype:** IgG  
**Applications:** ELISA, multiplex-based assays, and Western blot (WB); the recommended starting concentration for ELISA is 0.2-1 µg/ml, 0.5-2 µg/ml for WB, and 0.1-0.5 µg/ml for multiplex-based assays. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: Recombinant histone H3.3  
Lane 2: HeLa cells

**WB of Recombinant Histone H3.3 and Acid Extracts of HeLa Cells using 0.5 µg/ml of Histone H3K36Me2 Monoclonal Antibody (RM141).** This showed a band of histone H3 dimethylated at Lysine 36 (K36Me2) in HeLa cells.



Histone H3K36Me2 Monoclonal Antibody (RM141) Specifically Reacts to Histone H3 Dimethylates at Lysine 36 (K36Me2). No cross reactivity with unmodified H3K36, H3K36Me1, H3K36Me3, or other methylations in histone H3.

**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, 01/26/2024

**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 H3 is a nuclear protein and a component of the nucleosome core, a basic unit of chromatin, that is essential for organizing genomic DNA in eukaryotic nuclei.<sup>1</sup> It is a globular protein that contains an unstructured N-terminal tail that extends outside of the nucleosome core and is subject to various post-translational modifications (PTMs), including methylation, phosphorylation, acetylation, and citrullination.<sup>1,2</sup> Dimethylation of H3K36 is catalyzed by SET domain-containing histone methyltransferases including SET2, HYPB, NSD1, and ASH1L, and mutation of these methyltransferases is associated with various diseases, including multiple myeloma and Sotos syndrome.<sup>3-5</sup> H3K36Me2 is enriched at the coding regions of genes and correlates with the initiation, but not maintenance, of active transcription, heterochromatin maintenance, and DNA break repair.<sup>4,6,7</sup> Cayman's Histone H3K36Me2 Monoclonal Antibody (RM141) can be used for ELISA, multiplex-based assays, and Western blot (WB) applications.

## References

---

1. Hyun, K., Jeon, J., Park, K., *et al.* Writing, erasing and reading histone lysine methylations. *Exp. Mol. Med.* **49(4)**, e324 (2017).
2. Sharda, A., Amnekar, R.V., Natu, A., *et al.* Histone posttranslational modifications: Potential role in diagnosis, prognosis, and therapeutics of cancer. *Prognostic Epigenetics*. Sharma, S., editor, *Academic Press* (2019).
3. Li, J., Ahn, J.H., and Wang, G.G. Understanding histone H3 lysine 36 methylation and its deregulation in disease. *Cell Mol. Life Sci.* **76(15)**, 2899-2916 (2019).
4. Rao, B., Shibata, Y., Strahl, B.D., *et al.* Dimethylation of histone H3 at lysine 36 demarcates regulatory and nonregulatory chromatin genome-wide. *Mol. Cell. Biol.* **25(21)**, 9447-9459 (2005).
5. Sun, X.J., Wei, J., Wu, X.Y., *et al.* Identification and characterization of a novel human histone H3 lysine 36-specific methyltransferase. *J. Biol. Chem.* **280(42)**, 35261-35271 (2005).
6. Li, B., Jackson, J., Simon, M.D., *et al.* Histone H3 lysine 36 dimethylation (H3K36me2) is sufficient to recruit the Rpd3s histone deacetylase complex and to repress spurious transcription. *J. Biol. Chem.* **284(12)**, 7970-7976 (2009).
7. Avdic, V., Zhang, P., Lanouette, S., *et al.* Structural and biochemical insights into MLL1 core complex assembly. *Structure* **19(1)**, 101-108 (2011).

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