

PRODUCT INFORMATION

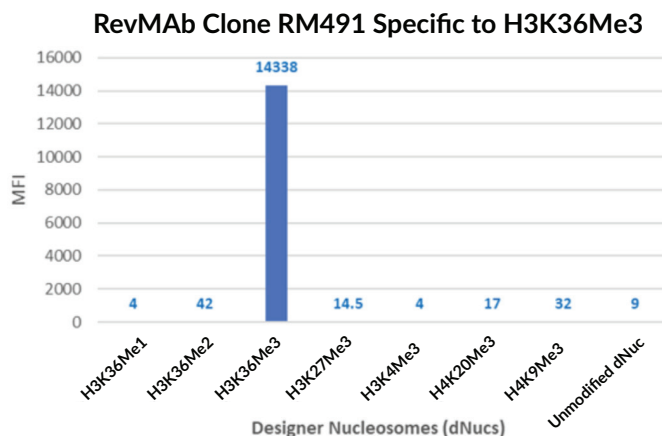
Histone H3K36Me3 Rabbit Monoclonal Antibody (Clone RM491)

Item No. 42795

Overview and Properties

Contents:	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
Synonym:	Trimethylated Histone 3 Lysine 36
Immunogen:	Peptide corresponding to H3K36Me3
Cross Reactivity:	(+) H3K36Me3; (-) H3K36Me1, H3K36Me2, H3K27Me3, H4K20Me3, H4K9Me3
Species Reactivity:	(+) Human, 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:	RM491
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA, Chromatin immunoprecipitation (ChIP), and multiplex-based assay; the recommended starting concentration is 0.2-1 µg/ml for ELISA, 2-10 µg/ml for ChIP, and 0.1-0.5 µg/ml for multiplex-based assay. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Luminex of Designer Recombinant Nucleosomes (dNucs) with Histone H3K36Me3 Rabbit Monoclonal Antibody (Clone RM491). Unmodified or modified dNucs containing lysine methylations were used.

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, 01/13/2025

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 3 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.¹ It is a globular protein that contains a histone fold domain with a C-terminal α -helix that facilitates nucleosome interactions and chromatin compaction, as well as an unstructured N-terminal tail that extends outside of the nucleosome core, both of which are subject to various post-translational modifications (PTMs), including ubiquitination, acetylation, methylation, and phosphorylation.¹⁻³ Histone H3 has a key role in transcriptional regulation.⁴ Increased nucleosome levels of trimethylated lysine 36 on histone 3 (H3K36Me3) are required for efficient DNA damage and nucleotide mismatch repair, DNA transcription, and repressing heterochromatin formation.⁵⁻⁷ Increased tumor levels of histone H3K36Me3 are correlated with increased tumor size, frequency of lymph node metastasis, and survival in patients with esophageal squamous cell carcinoma.⁸ Cayman's Histone H3K36Me3 Rabbit Monoclonal Antibody (Clone RM491) can be used for ELISA, chromatin immunoprecipitation (ChIP), and multiplex-based assay 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. Wyrick, J.J. and Parra, M.A. The role of histone H2A and H2B post-translational modifications in transcription: A genomic perspective. *Biochim. Biophys. Acta* **1789**(1), 37-44 (2009).
3. Wang, C.-Y., Hua, C.-Y., Hsu, H.-E., *et al.* The C-terminus of histone H2B is involved in chromatin compaction specifically at telomeres, independently of its monoubiquitylation at lysine 123. *PLoS One* **6**(7), e22209 (2011).
4. Bhaumik, S.R., Smith, E., and Shilatifard, A. Covalent modifications of histones during development and disease pathogenesis. *Nat. Struct. Mol. Biol.* **14**(11), 1008-1016 (2007).
5. Sun, Z., Zhang, Y., Jia, J., *et al.* H3K36me3, message from chromatin to DNA damage repair. *Cell Biosci.* **10**, 9 (2020).
6. Lucio-Eterovic, A.K., Singh, M.M., Gardner, J.E., *et al.* Role for the nuclear receptor-binding SET domain protein 1 (NSD1) methyltransferase in coordinating lysine 36 methylation at histone 3 with RNA polymerase II function. *Proc. Natl. Acad. Sci. USA* **107**(39), 16952-16957 (2010).
7. Harutyunyan, A.S., Chen, H., Lu, T., *et al.* H3K27M in gliomas causes a one-step decrease in H3K27 methylation and reduced spreading within the constraints of H3K36 methylation. *Cell Rep.* **33**(7), 108390 (2020).
8. Zhou, M., Li, Y., Lin, S., *et al.* H3K9me3, H3K36me3, and H4K20me3 expression correlates with patient outcome in esophageal squamous cell carcinoma as epigenetic markers. *Dig. Dis. Sci.* **64**(8), 2147-2157 (2019).

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