

# PRODUCT INFORMATION



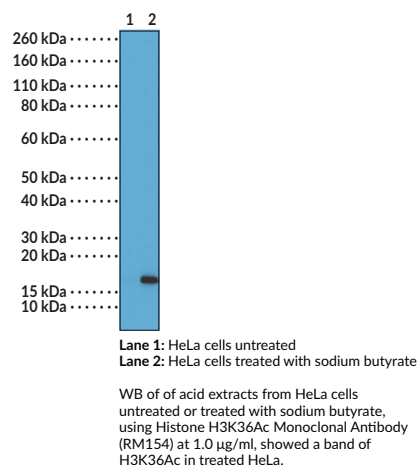
## Histone H3K36Ac Monoclonal Antibody (RM154)

Item No. 32135

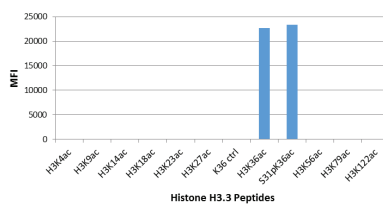
### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
<b>Synonyms:</b>	Acetylated Histone H3 Lysine 36
<b>Immunogen:</b>	Peptide corresponding to H3K36Ac
<b>Cross Reactivity:</b>	(+) H3K36Ac, H3K36Ac/S31Ph; (-) Unmodified H3K36, H3K4Ac, K3K9Ac, K3K14Ac, H3K18Ac, H3K23Ac, H3K27Ac, H3K56Ac, H3K79Ac, H3K122Ac
<b>Species Reactivity:</b>	(+) Vertebrates
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Concentration:</b>	1 mg/ml
<b>Clone:</b>	RM154
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	Chromatin immunoprecipitation (ChIP), ELISA, immunocytochemistry (ICC), multiplex-based assays, and Western blot (WB); the recommended starting concentration is 1-5 µg, 0.2-1 and 0.1-0.5 µg/ml for ChIP, ELISA, and multiplex-based assay, respectively, and 0.5-2 µg/ml for ICC and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

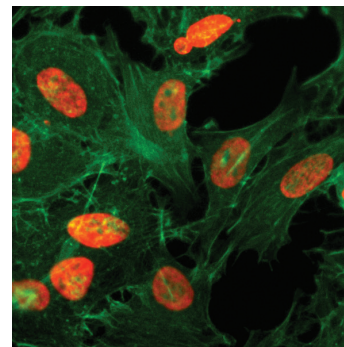
### Images



Histone H3K36Ac Monoclonal Antibody is Specific to Histone H3K36Ac



Histone H3K36Ac Monoclonal Antibody (RM 154) specifically reacts to histone H3 acetylated at lysine 36 (K36ac), and is not affected by the phosphorylation of neighboring Ser<sup>31</sup>. No cross reactivity with H3K4Ac, K3K9Ac, H3K14Ac, H3K18Ac, H3K23Ac, H3K27Ac, H3K56Ac, H3K79Ac, or H3K122Ac.



Immunocytochemistry of HeLa cells treated with sodium butyrate, using Histone H3K36Ac Monoclonal Antibody (RM154) (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, 02/07/2024

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> Acetylation of histone H3 at lysine 36 (H3K36Ac) is predominantly localized to the promoters of genes transcribed by RNA polymerase II and is highly conserved in mammals.<sup>3</sup> Cayman's Histone H3K36Ac Monoclonal Antibody (RM154) can be used for chromatin immunoprecipitation (ChIP), ELISA, immunocytochemistry (ICC), multiplex-based assay, 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. Morris, S.A., Rao, B., Garcia, B.A., *et al.* Identification of histone H3 lysine 36 acetylation as a highly conserved histone modification. *J. Biol. Chem.* **282(10)**, 7632-7640 (2007).

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