

# PRODUCT INFORMATION



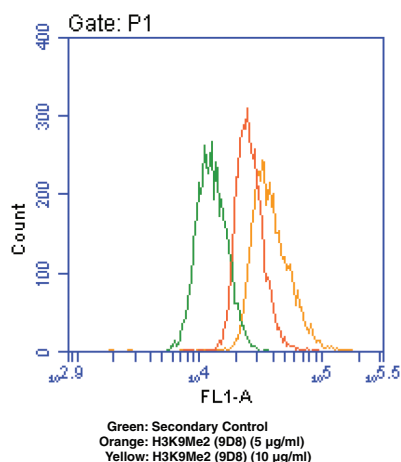
## Histone H3K9Me2 Monoclonal Antibody (Clone 9D8)

Item No. 18072

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of protein G-purified antibody.
<b>Synonym:</b>	Histone H3 containing dimethylated lysine 9
<b>Immunogen:</b>	Synthetic peptide from the N-terminal region of human H3K9Me2 conjugated to KLH
<b>Cross Reactivity:</b>	(+) H3K9Me3 (~50%); (-) H3K4Me2, H3K27Me2, H3K36Me2, H3K4Me
<b>Species Reactivity:</b>	(+) Human; other species not tested
<b>Uniprot No.:</b>	P68431
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol, 0.1% BSA, and 0.02% sodium azide
<b>Clone:</b>	9D8
<b>Host:</b>	Mouse
<b>Isotype:</b>	IgG2bk
<b>Applications:</b>	ELISA and Flow cytometry (FC); the recommended starting dilution for ELISA is 1:1,000 and 1:200 for FC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



HeLa cells were permeabilized and fixed using 4% formaldehyde. Cells were blocked with 1% FBS and probed with the indicated antibody.

**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, 11/20/2023

**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 histone H3 at lysine 9 (H3K9Me2) is associated with transcriptional repression.<sup>3</sup> Heterochromatin protein 1- $\alpha$  (HP1- $\alpha$ ), HP1- $\beta$ , and HP1- $\gamma$  selectively bind histone H3 N-terminal peptides containing dimethylated lysine 9 over unmodified lysine 9.<sup>4</sup> High nuclear levels of H3K9Me2 in tumor tissue are associated with decreased disease-specific and disease-free survival in patients with oral and oropharyngeal squamous cell carcinoma.<sup>5</sup> Cayman's Histone H3K9Me2 Monoclonal Antibody (Clone 9D8) can be used for ELISA and flow cytometry (FC) 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. Lienert, F., Mohn, F., Tiwari, V.K., *et al.* Genomic prevalence of heterochromatic H3K9me2 and transcription do not discriminate pluripotent from terminally differentiated cells. *PLoS Genet.* **7(6)**, e1002090 (2011).
4. Lachner, M., O'Carroll, D., Rea, S., *et al.* Methylation of histone H3 lysine 9 creates a binding site for HP1 proteins. *Nature* **410(6824)**, 116-120 (2001).
5. Maia, L.L., Peterle, G.T., Dos Santos, M., *et al.* JMJD1A, H3K9me1, H3K9me2 and ADM expression as prognostic markers in oral and oropharyngeal squamous cell carcinoma. *PLoS One* **13(3)**, e019884 (2018).

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