

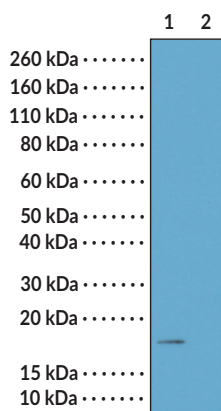
PRODUCT INFORMATION

Histone H3.3^{G34R} Rabbit Monoclonal Antibody - Biotinylated (Clone RM240)
Item No. 42868

Overview and Properties

Contents:	This vial contains 50 µl of protein A-affinity purified antibody.
Synonyms:	H3.3A, H3F3
Immunogen:	A peptide corresponding to histone H3.3 ^{G34R}
Cross Reactivity:	(+) Histone H3.3 ^{G34R} ; (-) Wild-type histone H3.3
Species Reactivity:	(+) Species Independent
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM240
Host:	Rabbit
Isotype:	IgG
Applications:	Chromatin Immunoprecipitation (ChIP), ELISA, Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1:250-1:2,500 for ELISA, 1:100-1:500 for IHC, an 1:200-1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: H3.3 G34R
Lane 2: H3.3 WT

WB of 293T cell lysates transfected with a DNA construct encoding wild-type or G34R mutant histone H3.3 using Histone H3.3^{G34R} Rabbit Monoclonal Antibody - Biotinylated (Clone RM240).

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, 03/12/2025

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 H3 is a nuclear protein and 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 an unstructured N-terminal tail that extends outside of the nucleosome core and is subject to various post-translational modifications (PTMs). Metazoans have three histone variants: canonical histone H3.1, DNA replication-independent histone H3.3, and the centromeric variant CENP-A.² Histone H3.3 is incorporated into chromatin in a DNA replication-independent manner at sites of active transcription.²⁻⁴ A missense mutation replacing guanine with thymine in *H3F3A*, the gene encoding histone H3.3, leads to a glycine-to-arginine substitution, histone H3.3^{G34R}, which is associated with glioblastoma multiforme in children and young adults.⁵ Cayman's Histone H3.3^{G34R} Rabbit Monoclonal Antibody (Clone RM240) - Biotinylated can be used for chromatin immunoprecipitation (ChIP), ELISA, immunohistochemistry (IHC), 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. Elsaesser, S.J., Goldberg, A.D., and Allis, C.D. New functions for an old variant: No substitute for histone H3.3. *Curr. Opin. Genet. Dev.* **20(2)**, 110-117 (2010).
3. Bjerke, L., Mackay, A., Nandhabalan, M., *et al.* Histone H3.3 mutations drive pediatric glioblastoma through upregulation of MYCN. *Cancer Discov.* **3(5)**, 512-519 (2013).
4. Tagami, H., Ray-Gallett, D., Almouzni, G., *et al.* Histone H3.1 and H3.3 complexes mediate nucleosome assembly pathways dependent or independent of DNA synthesis. *Cell* **116(1)**, 51-61 (2004).
5. Schwartzenuber, J., Korshunov, A., Liu, X.-Y., *et al.* Driver mutations in histone H3.3 and chromatin remodelling genes in paediatric glioblastoma. *Nature* **482(7384)**, 226-231 (2012).

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