

PRODUCT INFORMATION



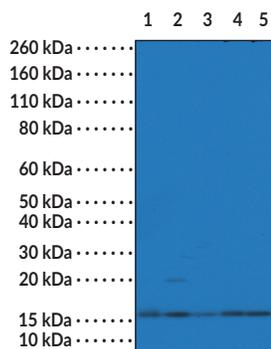
Histone H2A.Z (C-Term) Monoclonal Antibody (Clone RM215)

Item No. 32184

Overview and Properties

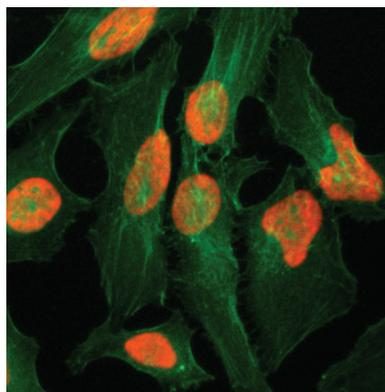
Contents:	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
Synonym:	H2AZ
Immunogen:	Peptide from the C-terminal region of human histone H2A.Z
Cross Reactivity:	(+) H2A.Z independent of PTMs; (-) Other histone proteins
Species Reactivity:	(+) Vertebrates
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Concentration:	1.0 mg/ml
Clone:	RM215
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA, Immunocytochemistry (ICC), Multiplex-based assay, and Western blot (WB); the recommended starting concentration is 0.1-1 µg/ml for ELISA and multiplex-based assays, 1-2 µg/ml for ICC, and 0.1-0.5 µg/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: HeLa cell lysates
Lane 2: HEK293 cell lysates
Lane 3: A375 cell lysates
Lane 4: SK-MEL-2 cell lysates
Lane 5: A431 cell lysates

WB of A375, A431, HEK293, HeLa, and SK-MEL-2 whole cell lysates, using Histone H2A.Z (C-Term) Monoclonal Antibody (Clone RM215) at a concentration of 0.5 µg/ml.



Immunofluorescent labeling of HeLa cells, using Histone H2A.Z (C-Term) Monoclonal Antibody (Clone RM215) (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, 01/18/2024

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 H2A.Z is a variant of histone H2A, a nuclear protein and a component of the nucleosome core.^{1,2} It is a globular protein containing unstructured N- and C-terminal tails that extend outside of the nucleosome core that are subject to a variety of post-translational modifications (PTMs), including acetylation, methylation, and ubiquitination, which function as epigenetic regulators of transcription.³ Histone H2A.Z influences chromatin accessibility and nucleosome stability, regulating various cellular functions, including DNA replication and repair, mitosis, heterochromatin formation, and chromosome segregation.^{1,4} It is enriched at gene promoters and associated with active or repressed gene transcription in a manner dependent on species or PTMs, among other factors.^{4,5} Increased tumor H2A.Z levels are associated with decreased overall survival in patients with breast cancer.⁶ Cayman's Histone H2A.Z (C-Term) Monoclonal Antibody (Clone RM215) can be used for ELISA, immunocytochemistry (ICC), multiplex-based assay, and Western blot (WB) applications. The antibody recognizes the C-terminal region of histone H2A.Z independent of PTMs.

References

1. Bönisch, C. and Hake, S.B. Histone H2A variants in nucleosomes and chromatin: More or less stable? *Nucleic Acids Res.* **40(21)**, 10719-10741 (2012).
2. Hyun, K., Jeon, J., Park, K., *et al.* Writing, erasing and reading histone lysine methylations. *Exp. Mol. Med.* **49(4)**, e324 (2017).
3. Corujo, D. and Buschbeck, M. Post-translational modifications of H2A histone variants and their role in cancer. *Cancers (Basel)* **10(3)**, 59 (2018).
4. Monteiro, F.L., Baptista, T., Amado, F., *et al.* Expression and functionality of histone H2A variants in cancer. *Oncotarget* **5(11)**, 3428-3443 (2014).
5. Marques, M., Laflamme, L., Gervais, A.L., *et al.* Reconciling the positive and negative roles of histone H2A.Z in gene transcription. *Epigenetics* **5(4)**, 267-272 (2010).
6. Hua, S., Kallen, C.B., Dhar, R., *et al.* Genomic analysis of estrogen cascade reveals histone variant H2A.Z associated with breast cancer progression. *Mol. Syst. Biol.* **4(1)**, 188 (2008).

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