

PRODUCT INFORMATION



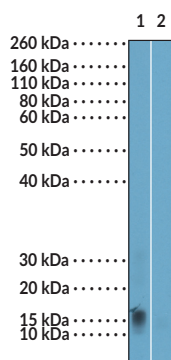
Histone H2BK11Ac Rabbit Monoclonal Antibody (Clone RM456)

Item No. 42791

Overview and Properties

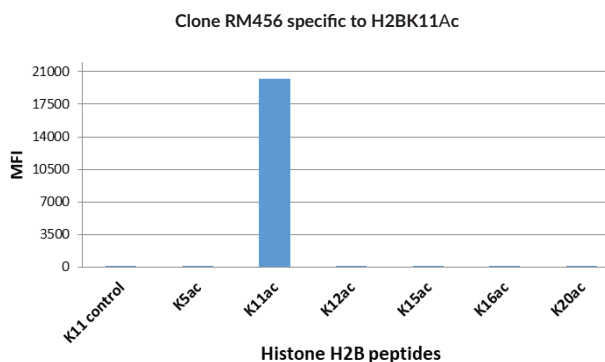
Contents:	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
Synonyms:	Acetylated Histone H2B Lysine 11
Immunogen:	An acetyl-peptide corresponding to Acetyl-Histone H2B (Lys11)
Cross Reactivity:	(+) H2BK11Ac; (-) Unmodified H2BK11, H2B5Ac, H2BK12Ac, H2BK15Ac, H2BK16Ac, H2BK20Ac
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
Clone:	RM456
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA, multiplex-based assay, and Western blot (WB); the recommended starting concentration is 0.2-1 µg/ml for ELISA, 0.01-1 µg/ml for multiplex-based assay, and 0.001-0.01 µg/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: HeLa cells treated
Lane 2: HeLa cells untreated

WB of HeLa cells treated with sodium butyrate or left untreated using Histone H2BK11Ac Rabbit Monoclonal Antibody (Clone RM456) at a concentration of 0.002 µg/ml.



Histone H2BK11Ac Rabbit Monoclonal Antibody (Clone RM456) specifically reacts to histone H2B acetylated at Lysine 11 (K11Ac). No cross reactivity with unmodified Lysine 11, or other acethylations in histone H2B.

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, 01/10/2025

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 H2B 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 Acetylation of histone H2B at lysine 11 (H2BK11Ac) is downregulated in a mouse model of allergic asthma.⁴ Cayman's Histone H2BK11Ac Rabbit Monoclonal Antibody (Clone RM456) can be used for ELISA, 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. 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. Ren, Y., Li, M., Bai, S., *et al.* Identification of histone acetylation in a murine model of allergic asthma by proteomic analysis. *Exp. Biol. Med. (Maywood)* **246**(8), 929-939 (2021).

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