

PRODUCT INFORMATION



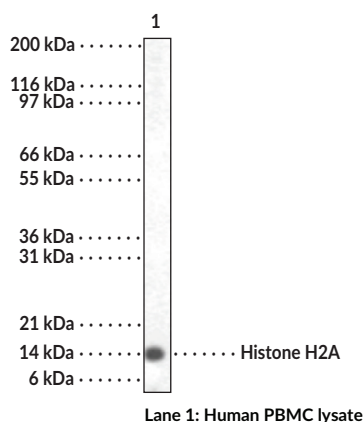
Histone H2A Polyclonal Antibody

Item No. 13535

Overview and Properties

Contents:	100 µg of protein G-purified IgG in 200 µl PBS with 0.05% BSA and 0.05% sodium azide
Immunogen:	Synthetic peptide that corresponds to amino acids 1-15 (MSGRGKQGKARAKA) and 81-96 (PRHLQLAIRNDEELNK) in human histone H2A
Cross Reactivity:	(+) Histone H2A
Species Reactivity:	(+) Human, mouse
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Host:	Rabbit
Applications:	ELISA and Western blot (WB); the recommended starting dilution is 1:1,000 for ELISA and 1 µg/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



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, 01/31/2022

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 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 an unstructured N-terminal tail that extends outside of the nucleosome core and is subject to a variety of post-translational modifications (PTMs), including ubiquitination, acetylation, methylation, and phosphorylation, which function as epigenetic regulators of transcription.^{1,2} Serum histone H2A autoantibodies have been found in patients with drug-induced or spontaneous systemic lupus erythematosus (SLE).³ Cayman's Histone H2A Polyclonal Antibody can be used for Western blot (WB) and ELISA 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. Wang, Y., Li, M., Stadler, S., *et al.* Histone hypercitrullination mediates chromatin decondensation and neutrophil extracellular trap formation. *J. Cell Biol.* **184(2)**, 205-213 (2009).
3. Burlingame, R.W. and Rubin, R.L. Autoantibody to the nucleosome subunit (H2A-H2B)-DNA is an early and ubiquitous feature of lupus-like conditions. *Mol. Biol. Rep.* **23(3-4)**, 159-166 (1996).

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