

PRODUCT INFORMATION



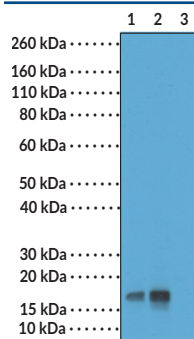
Histone H3T3Ph Monoclonal Antibody (RM159)

Item No. 32165

Overview and Properties

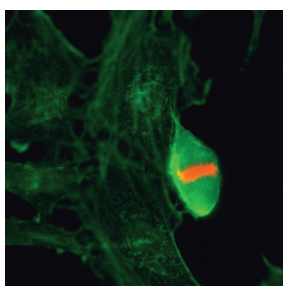
Contents:	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
Synonyms:	H3pT3, H3pThr3, Histone H3 (Phospho-Thr3), Phospho-Histone H3 Threonine 3, Phosphorylated Histone H3 Threonine 3
Immunogen:	Peptide corresponding to H3T3Ph
Cross Reactivity:	(+) Vertebrates
Species Reactivity:	(+) H3T3Ph; (-) Other phosphorylated histones
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:	RM159
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA, immunocytochemistry (ICC), immunohistochemistry (IHC), multiplex-based assays, and Western blot (WB); the recommended starting concentration is 0.2-1 µg/ml for ELISA, 0.5-2 µg/ml for ICC and IHC, and 0.1-1 µg/ml for multiplex-based assays and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images

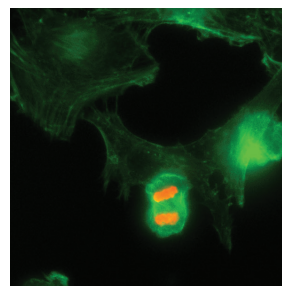


Lane 1: Acid extracts of HeLa cells untreated
Lane 2: Acid extracts of HeLa cells treated
Lane 3: Recombinant histone H3.3

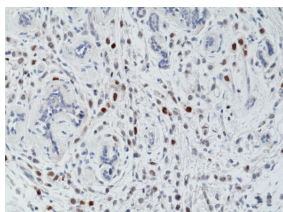
WB of acid extracts of HeLa cells treated with nocodazole or left untreated, and recombinant histone H3.3 protein using Histone H3T3Ph Monoclonal Antibody (RM159) at a concentration of 0.1 µg/ml.



Immunocytochemical staining of HeLa cells using Histone H3T3Ph Monoclonal Antibody (RM159) (red). Actin filaments have been labeled with fluorescein phalloidin (green).



Immunocytochemical staining of HeLa cells using Histone H3T3Ph Monoclonal Antibody (RM159) (red). Actin filaments have been labeled with fluorescein phalloidin (green).



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast cancer tissue using Histone H3T3Ph Monoclonal Antibody (RM159) at a concentration of 1 µg/ml.

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, 02/29/2024

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 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.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), including methylation, phosphorylation, acetylation, and citrullination.^{1,2} Phosphorylation of histone H3 at threonine 3 (H3T3Ph), catalyzed by the serine/threonine protein kinase haspin, is involved in the localization of the chromosomal passenger complex (CPC) at centromeres during mitosis and chromatin condensation during meiosis.^{3,4} Cayman's Histone H3T3Ph Monoclonal Antibody (RM159) can be used for ELISA, immunocytochemistry (ICC), immunohistochemistry (IHC), 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. 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. Kelly, A.E., Ghenoiu, C., Xue, J.Z., *et al.* Survivin reads phosphorylated histone H3 threonine 3 to activate the mitotic kinase Aurora B. *Science* **330(6001)**, 235-239 (2010).
4. Wang, Q., Wei, H., Du, J., *et al.* H3 Thr3 phosphorylation is crucial for meiotic resumption and anaphase onset in oocyte meiosis. *Cell Cycle* **15(2)**, 213-224 (2016).

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