

PRODUCT INFORMATION

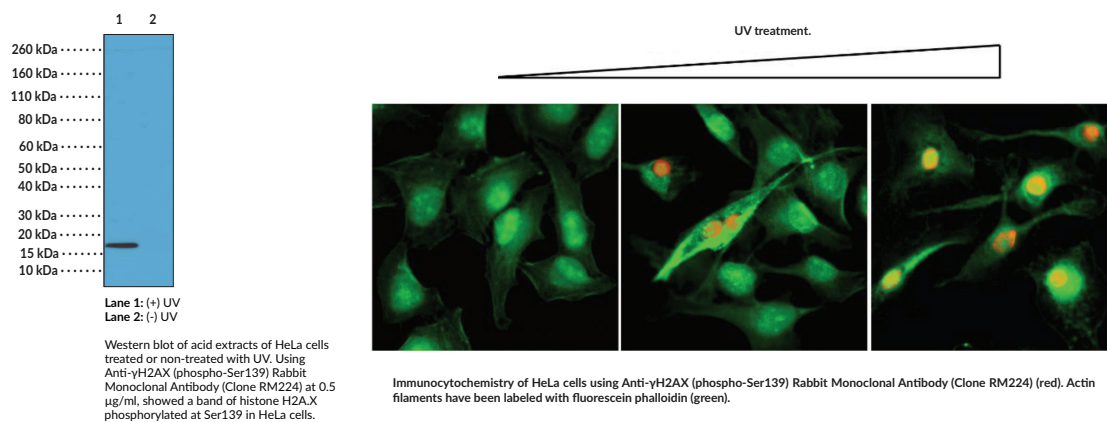
Anti- γ H2AX (phospho-Ser139) Rabbit Monoclonal Antibody (Clone RM224)

Item No. 20721

Overview and Properties

Contents:	This vial contains 100 μ g of protein A affinity-purified antibody from an animal origin-free culture supernatant.
Immunogen:	A phospho-peptide corresponding to γ H2AX (phospho-Ser139)
Cross Reactivity:	(+) Phospho- γ H2AXSer139; (-) Other phosphorylated histones
Species Reactivity:	(+) All species
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	\geq 1 year
Storage Buffer:	50% glycerol/PBS with 1% BSA and 0.09% sodium azide
Clone:	RM224
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA, Immunocytochemistry (ICC), Multiplex, and Western blot (WB); the recommended starting dilution for ELISA is 0.2 - 1 μ g/ml, 0.5 - 2 μ g/ml for ICC and WB, and 0.1 - 1 μ g/ml for multiplex. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



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, 01/10/2022

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 H2AX, a variant of the nucleosome core histone H2A, undergoes phosphorylation on serine 139 in response to DNA damage, particularly when the damage involves formation of double-strand breaks. Phosphorylated H2AX is also termed γ H2AX. Induction of γ H2AX in cells exposed to genotoxic agents is considered a sensitive and specific reporter of DNA damage.¹ This antibody reacts to γ H2AX when phosphorylated at Ser139 and does not detect other phosphorylated histones.

Reference

1. Huang, X., Tanaka, T., Kurose, A., *et al.* Constitutive histone H2AX phosphorylation on Ser-139 in cells untreated by genotoxic agents is cell-cycle phase specific and attenuated by scavenging reactive oxygen species. *Int. J. Oncol.* **29(2)**, 495-501 (2006).

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