

[pSer¹³⁹]Histone H2AX monoclonal antibody (9F3)

The nucleosome is made up of four core histone proteins (H2A, H2B, H3 and H4) and is the primary building block of chromatin. The core histones form an octamer, which is made up of one H3-H4 tetramer and two H2A-H2B dimers. The N-terminal tail of core histones undergoes different post-transcriptional modifications including acetylation, phosphorylation, methylation and ubiquitination. These modifications occur in response to cell signal stimuli and have a direct effect on gene expression.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

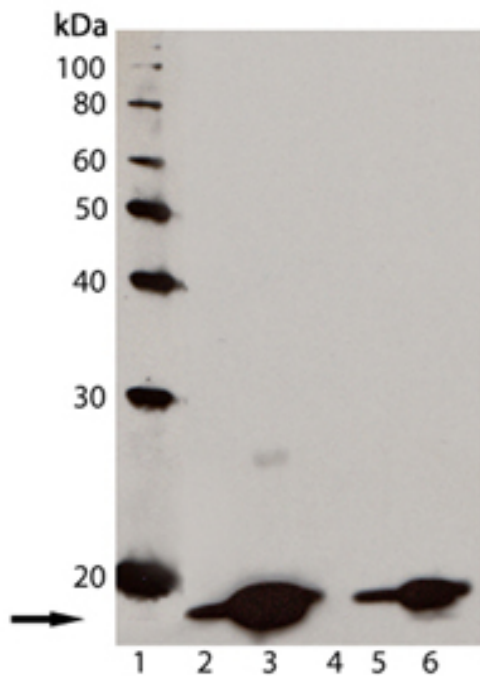
Ordering Information

[Order Online »](#)

ADI-KAM-CC255-D	50µg
ADI-KAM-CC255-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of [pSer139] Histone H2AX, mAb (9F3) (Prod. No. ADI-KAM-CC255): Lane 1: MW marker; Lane 2: Jurkat cell lysate; Lane 3: Jurkat cell lysate treated with staurosporine; Lane 4: 3T3 cell lysate; Lane 5: CHO-K1 cell lysate; Lane 6: Rat-2 cell lysate.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

<h2>Product Details</h2>	
Application	WB
Application Notes	Detects a band of ~16kDa by Western blot.
Clone	9F3
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to a portion of human Histone H2AX phosphorylated at Ser139.
Isotype	IgG1k
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites.
Species Reactivity	Bovine, Chicken, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep
Technical Info / Product Notes	Patents: USPN 6,362,317 and 6,884,873
UniProt ID	P16104

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)