

# [pSer<sup>139</sup>]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 »](#)

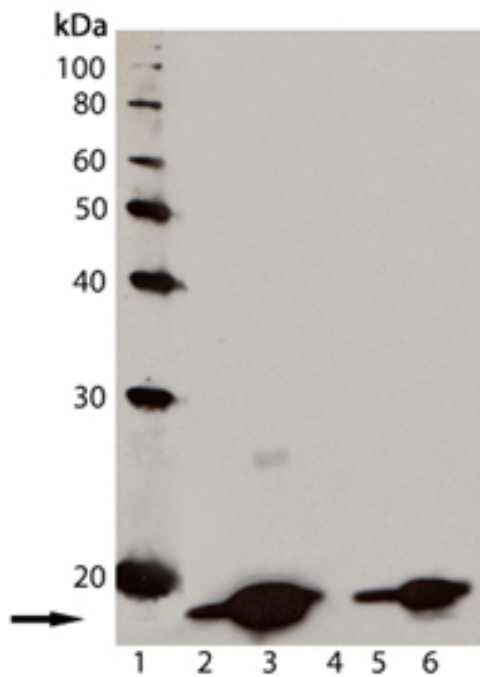
## Ordering Information

[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

## Product Details

**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** IgG1κ

**Purity Detail** Protein G affinity purified.

**Recommendation** 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**



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  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

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

**Worry-free Guarantee**

**This antibody is covered by our Worry-Free Guarantee.**