# [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.

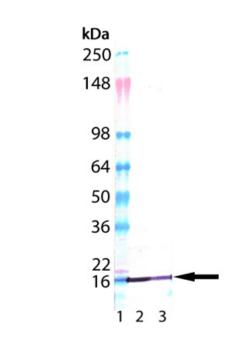
## **Ordering Information**

**Order Online** »

ENZ-ABS662-0050	50µg
ENZ-ABS662-0200	200μg

Manuals, SDS & CofA

**View Online** »



Western Blot analysis of ENZ-ABS662: Lane 1: MW Marker, Lane 2: Jurkat cell lysate treated with staurosporine; Lane 3: 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

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 IgG1ĸ

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 hybridoma tissue culture supernatant.

**Species Reactivity**Bovine, Chicken, Dog, Guinea pig, Hamster, Human,

Monkey, Mouse, Porcine, Rabbit, Rat

UniProt ID P16104

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com