[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 »

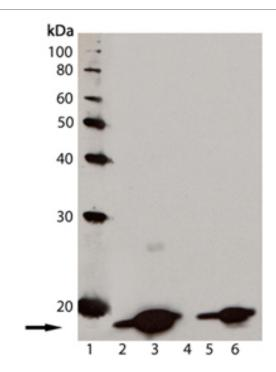
Ordering Information

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

Product Details

Application WB

Application NotesDetects 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.

lgG1k

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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com