

Histone H2A antibody (pAb)

Catalog Nos: 39235, 39436, 39236

RRID: AB_2687477

Isotype: Serum

Application(s): ChIP, IP, WB

Reactivity: Budding Yeast, Fission Yeast

Volumes: 100 µl, 50 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 14 kDa

Background: Histone H2A is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 147 base pairs of DNA wrapped around an octamer of core histone proteins (two each of Histone H2A, Histone H2B, Histone H3 and Histone H4). Histone H1 is a linker histone, present at the interface between the nucleosome core and DNA entry/exit points; it is responsible for establishing higher-order chromatin structure. Chromatin is subject to a variety of chemical modifications, including post-translational modifications of the histone proteins and the methylation of cytosine residues in the DNA. Reported histone modifications include acetylation, methylation, phosphorylation, ubiquitylation, glycosylation, ADP-ribosylation, carbonylation and SUMOylation; they play a major role in regulating gene expression.

Immunogen: This Histone H2A antibody was raised against recombinant yeast histone H2A.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic. For your convenience, an IgG version (Catalog No. 39945) of this antibody that was purified by Protein A Chromatography is also available.

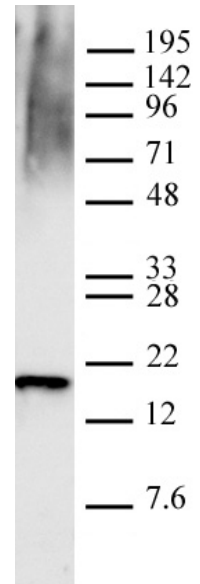
Application Notes:

Applications Validated by Active Motif:

WB: 1:2,000 - 1:5,000 dilution

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Histone H2A antibody (pAb) tested by Western blot.

Detection of Histone H2A by Western blot. The analysis was performed using *S. cerevisiae* whole cell acid extract probed with Histone H2A antibody at a 1:4,000 dilution.