

## Histone H2A antibody (pAb)

---

**Catalog No:** 39209

**RRID:** AB\_2793184

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Quantity:** 100 µg

**Purification:** Affinity Purified

**Host:** Rabbit

**Concentration:** 1 µg/µl

**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 a mixture of synthetic peptides corresponding to amino acid residues 1-5 and 81-96 of human Histone H2A.

**Buffer:** PBS containing 0.02% sodium azide and 0.05% BSA. Sodium azide is highly toxic.

**Application Notes:**

Validated Applications:

WB: 2 µg/ml dilution

Published Applications: see references

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all Blotto solutions may also reduce background. Individual optimization may be required.

**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 pAb tested by Western blot.**

Detection of Histone H2A by Western blot. The analysis was performed using PBMC whole-cell extract and Histone H2A pAb at a 2 µg/ml dilution.