

Histone H1 antibody (pAb)

Catalog No: 39575

Isotype: IgG

Application(s): ChIP, IF, IP, WB

Reactivity: Drosophila

Purification: Affinity Purified

Host: Rabbit

Volume: 100 µl

Molecular Weight: 38 kDa

Background: 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. Histone H1 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 and play a major role in regulating gene expression.

Immunogen: This Histone H1 antibody was raised against a recombinant protein corresponding to full-length *Drosophila* Histone H1.

Buffer: PBS containing 0.1% sodium azide, 0.2 mg/ml BSA and 35% glycerol. Sodium azide is highly toxic.

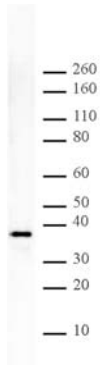
Application Notes: This Histone H1 antibody has been validated for use in Western blotting (1:2,000 - 1:10,000 dilution), immunofluorescence (1:1,000 - 1:2,500 dilution) and immunoprecipitation (5 µl per IP). Addition of 0.05% Tween 20 in blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot.

References: This antibody has been used in the following references:

ChIP: Lu X *et al.* (2009) *Genes Dev* 23: 452-465.

Storage and Guarantee: Antibodies in solution can be stored at -20°C for 2 years. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

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

**Histone H1 pAb tested by Western blot.**

Detection of Histone H1 by Western blot. The analysis was performed using Schneider's *Drosophila* L2 nuclear extract and Histone H1 pAb at a 1:5,000 dilution.