

Histone H1 antibody (mAb)

Catalog Nos: 61785, 61786

RRID: AB_2793765

Clone: 1G6-E3

Isotype: IgG1

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

Reactivity: Drosophila

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 35 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 antibody was raised against a recombinant protein containing amino acids 109-256 of *Drosophila melanogaster* Histone H1. This region of the protein is not conserved in mammals; the antibody is not likely to cross-react to human, mouse or rat Histone H1.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

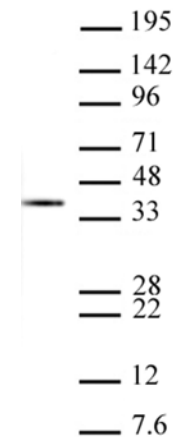
Application Notes:

Applications Validated by Active Motif:

WB: 0.2 - 1 µg/ml

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 H1 antibody tested by Western Blot.
Schneider's L2 *Drosophila* cell lysate (cat. no. 36087) probed with Histone H1 antibody (0.5 µg/ml)