

SUV39H1 antibody (mAb)

Catalog Nos: 39785, 39786

RRID: AB_2793343

Clone: MG44

Isotype: IgG1

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

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 50 kDa

Background: SUV39H1 is one of two homologues of the *Drosophila* Suppressor-of-variegation 3(9) protein (aka Su(var)3-9). SUV39H1 is a SET domain-containing histone methyltransferase specific for lysine 9 of histone H3. Lysine 9 methylation of histone H3 is found at regions of the genome that are transcriptionally silenced, such as centromeres, repetitive elements and inactive genes. H3 lysine 9 methylation is a hallmark of heterochromatin, regions of the genome that are in a tightly compacted state and relatively devoid of gene expression. HP1 proteins recognize and bind lysine 9 methylation and this contributes to maintaining heterochromatin in its compact and inactive state. SUV39H1 also contains a chromodomain, which may bind to lysine 9 methylation and contribute to the regional spreading of lysine 9 methylation and the establishment of heterochromatin over broader domains of the genome.

Immunogen: This SUV39H1 antibody was raised against a recombinant protein corresponding to amino acids 1-126 of mouse SUV39H1.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

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.