
SUV39H2 Polyclonal Antibody

(Catalog # A-3010)

Background

SUV39H2 encodes a H3 HMTase that shares 59% identity with SUV39H1, but which differs by the presence of a highly basic N terminus. SUV39H2 is also responsible for creating HP1 binding sites in heterochromatin by methylation of histone H3-Lys9. Moreover, SUV39H2 specifically accumulates with chromatin of the sex chromosomes (XYbody), which undergo transcriptional silencing during the first meiotic prophase.

Description

SUV39H2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q9H511

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 141-350 of human SUV39H2 (NP_001180354.1).

Storage

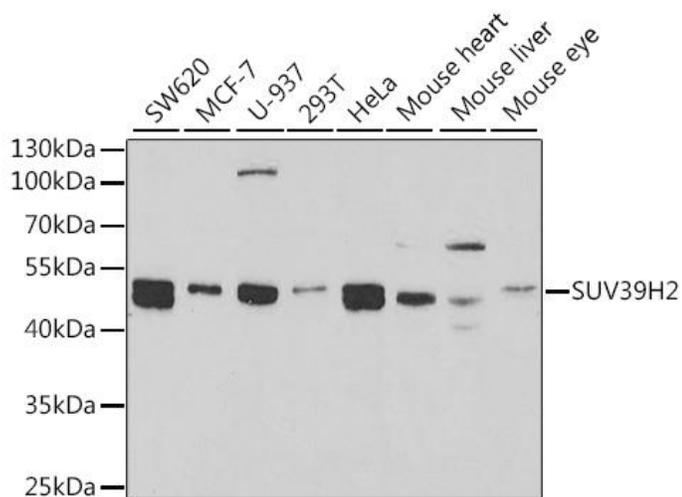
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

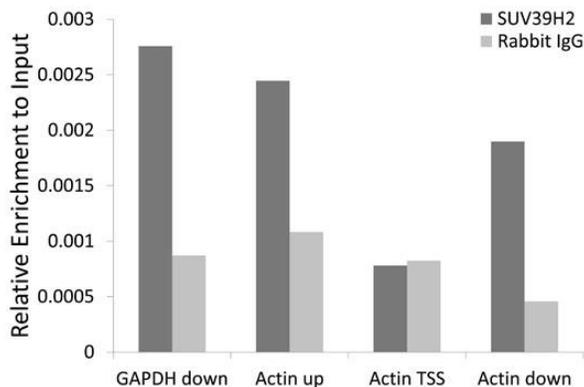
FLJ23414 antibody, H3 K9 HMTase 2 antibody, Histone H3 K9 methyltransferase 2 antibody, Histone lysine N methyltransferase H3 lysine 9 specific 2 antibody, Histone lysine N methyltransferase SUV39H2 antibody, KMT1B antibody, Lysine N methyltransferase 1B antibody, Suppressor of variegation 3 9 homolog 2 (Drosophila) antibody, Su(var)3 9 Drosophila homolog of 2 antibody, Suppressor of variegation 3 9 homolog 2 antibody, Suv39h2 antibody, SUV92_HUMAN antibody

Application

WB, IF, ChIP; Recommended dilution: WB 1:500 - 1:2000, IF 1:50 - 1:200, ChIP 1:20 - 1:100



Western blot analysis of extracts of various cell lines, using SUV39H2 antibody at 1:1000 dilution.
 Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.



Chromatin immunoprecipitation analysis of extracts of 293T cell line, using SUV39H2 antibody and rabbit IgG. The amount of immunoprecipitated DNA was checked by quantitative PCR. Histogram was constructed by the ratios of the immunoprecipitated DNA to the input.