
Histone H3K27me2 (H3K27 Dimethyl) Polyclonal Antibody

(Catalog # A-4038)

Background

Modulation of chromatin structure plays an important role in the regulation of transcription in eukaryotes. The nucleosome, made up of DNA wound around eight core histone proteins (two each of H2A, H2B, H3, and H4), is the primary building block of chromatin. The amino-terminal tails of core histones undergo various post-translational modifications, including acetylation, phosphorylation, methylation, and ubiquitination. These modifications occur in response to various stimuli and have a direct effect on the accessibility of chromatin to transcription factors and, therefore, gene expression. In most species, histone H2B is primarily acetylated at Lys5, 12, 15, and 20. Histone H3 is primarily acetylated at Lys9, 14, 18, 23, 27, and 56. Acetylation of H3 at Lys9 appears to have a dominant role in histone deposition and chromatin assembly in some organisms. Phosphorylation at Ser10, Ser28, and Thr11 of histone H3 is tightly correlated with chromosome condensation during both mitosis and meiosis. Phosphorylation at Thr3 of histone H3 is highly conserved among many species and is catalyzed by the kinase haspin. Immunostaining with phospho-specific antibodies in mammalian cells reveals mitotic phosphorylation at Thr3 of H3 in prophase and its dephosphorylation during anaphase.

Description

Histone H3K27me2 (H3K27 Dimethyl) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Broad Range, Mouse, Rat, Human

Isotype

IgG

Uniprot ID

Q16695

Purification

Affinity Purified

Immunogen

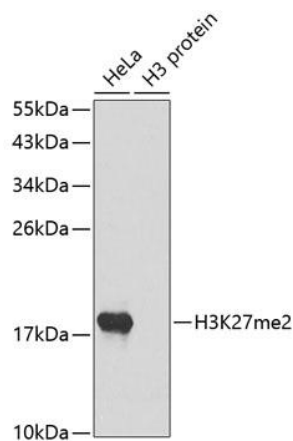
Synthetic Peptide of Human DiMethyl-Histone H3-K27

Storage

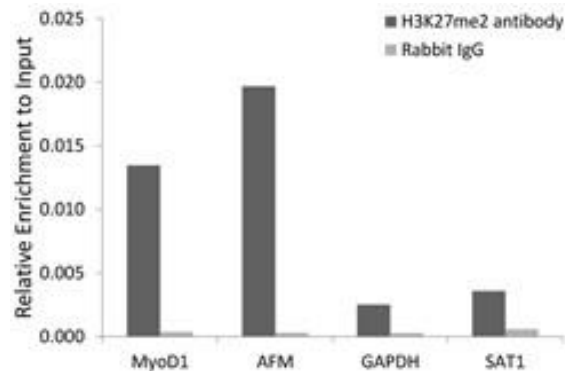
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze.

Application

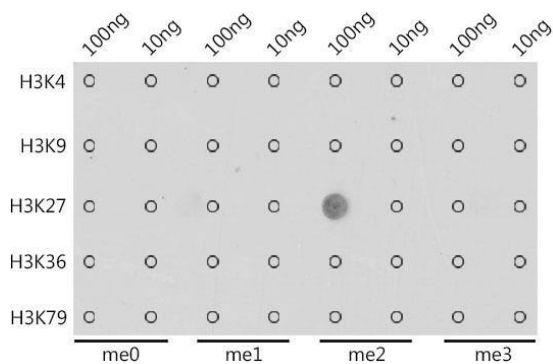
WB, IF, IP, ChIP, ChIP-seq; Recommended dilution: WB 1:500 - 1:2000, IF 1:50 - 1:200, IP 1:50 - 1:200, ChIP 1:20 - 1:100, ChIP-seq 1:20 - 1:100



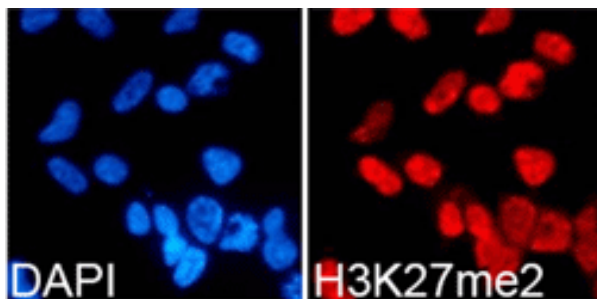
Western blot analysis of extracts of various cell lines, using DiMethyl-Histone H3-K27 antibody.
 Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.



Chromatin immunoprecipitation analysis extracts of 293 cell line, using Histone H3K27 Dimethyl Polyclonal 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.



Dot-blot analysis of all sorts of methylation peptides using DiMethyl-Histone H3-K27 antibody at 1:1000 dilution.



Immunofluorescence analysis of 293T cell using H3K27me2 Dimethyl Polyclonal Antibody. Blue: DAPI for nuclear staining.