

Histone H3K36me2 (H3K36 Dimethyl) Polyclonal Antibody

(Catalog # A-4041)

Background

Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. Nucleosomes consist of approximately 146 bp of DNA wrapped around a histone octamer composed of pairs of each of the four core histones (H2A, H2B, H3, and H4). The chromatin fiber is further compacted through the interaction of a linker histone, H1, with the DNA between the nucleosomes to form higher order chromatin structures. This gene is intronless and encodes a replication-dependent histone that is a member of the histone H3 family. Transcripts from this gene lack polyA tails; instead, they contain a palindromic termination element. This gene is located separately from the other H3 genes that are in the histone gene cluster on chromosome 6p22-p21.3.

Description

Histone H3K36me2 (H3K36 Dimethyl) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.05% proclin300 ,50% glycerol, pH7.3.

Specificity

Broad Range, Mouse, Rat, Human

Isotype

IgG

Uniprot ID

Q16695/P68431

Purification

Affinity Purified

Immunogen

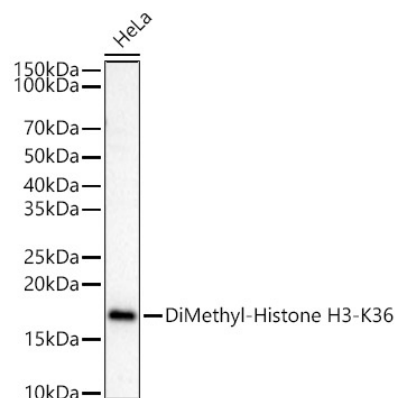
A synthetic dimethylated peptide around K36 of human histone H3 (NP_003520.1)

Storage

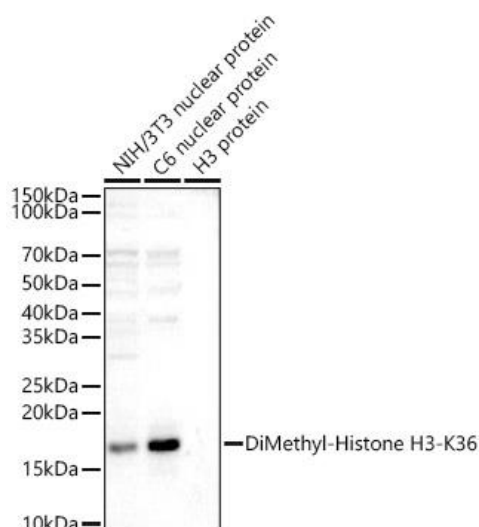
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Application

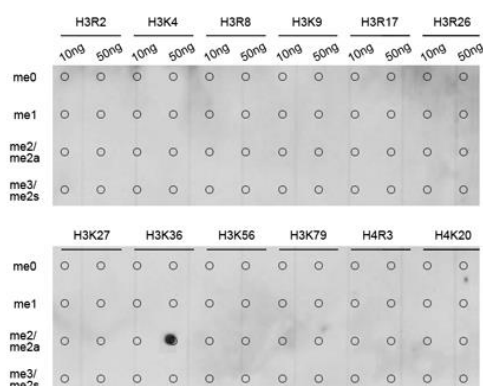
WB, IHC, IF/ICC, ChIP, ChIPseq, ELISA; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:200, IF/ICC 1:50 - 1:200, ChIP 5µg antibody for 5µg-10µg of Chromatin, CHIPseq 1:20 - 1:100, ELISA - recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements



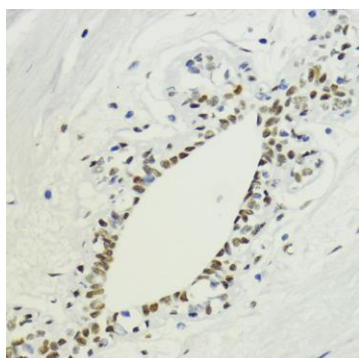
Western blot analysis of lysates from HeLa cells, using Histone H3K36me2 (H3K36 Dimethyl) Polyclonal Antibody at 1:600 dilution. Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25µg per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 180s.



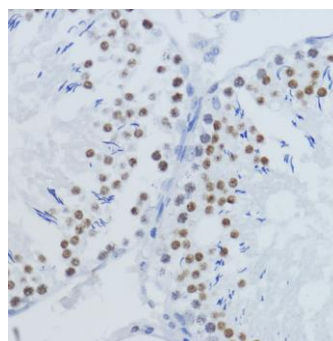
Western blot analysis of various lysates, using Histone H3K36me2 (H3K36 Dimethyl) Polyclonal Antibody at 1:600 dilution. Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25µg per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 180s.



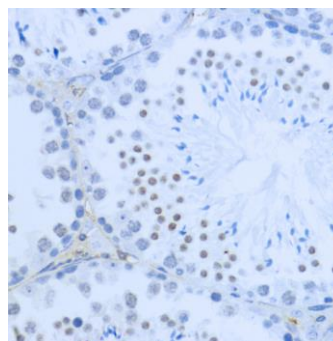
Dot-blot analysis of all sorts of methylation peptides using Histone H3K36me2 (H3K36 Dimethyl) Polyclonal Antibody.



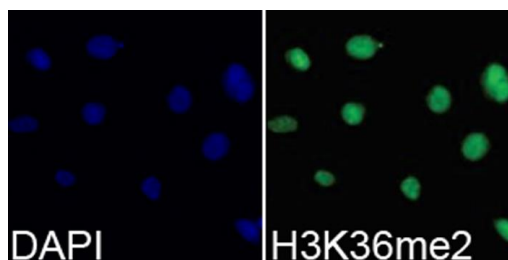
Immunohistochemistry of paraffin-embedded human breast using Histone H3K36me2 (H3K36 Dimethyl) Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



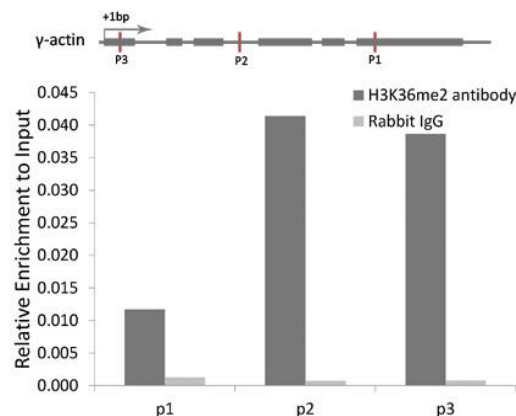
Immunohistochemistry of paraffin-embedded rat testis using Histone H3K36me2 (H3K36 Dimethyl) Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



Immunohistochemistry of paraffin-embedded mouse testis using Histone H3K36me2 (H3K36 Dimethyl) Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



Immunofluorescence analysis of 293T cells using Histone H3K36me2 (H3K36 Dimethyl) Polyclonal Antibody. Blue: DAPI for nuclear staining.



Chromatin immunoprecipitation analysis of γ -actin gene from 293 cell line, using Histone H3K36me2 (H3K36 Dimethyl) Polyclonal Antibody and rabbit IgG. P1, P2 and P3 were probes located on γ -actin gene as the schematic diagram illustrated. The amount of immunoprecipitated DNA was checked by quantitative PCR. Histogram was constructed by the ratios of the immunoprecipitated DNA to the input.