

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

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

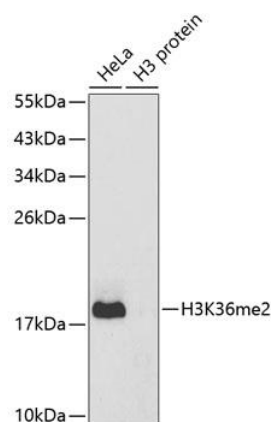
A synthetic dimethylated peptide around K36 of human histone H3 (NP_003520.1)

Storage

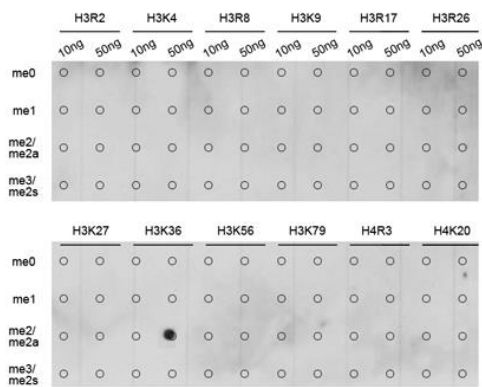
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Application

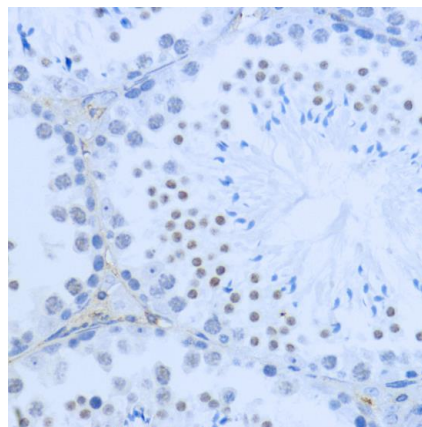
WB, IHC, IF, IP, ChIP, ChIPseq; Recommended dilution: WB 1:500 - 1:2000 IHC 1:50 - 1:200 IF 1:50 - 1:200 IP 1:50 - 1:200 ChIP 1:20 - 1:100 CHIPseq 1:20 - 1:100



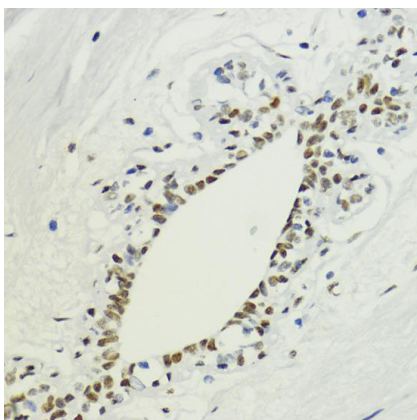
Western blot analysis of extracts of various cell lines, using DiMethyl-Histone H3-K36 antibody.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.



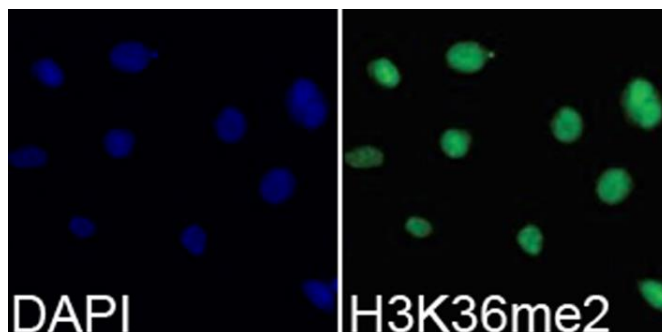
Dot-blot analysis of all sorts of methylation peptides using DiMethyl-Histone H3-K36 antibody.



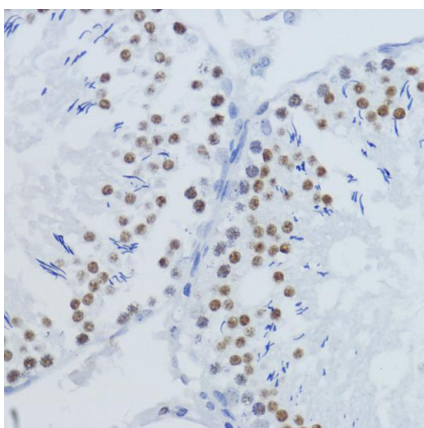
Immunohistochemistry of paraffin-embedded mouse testis using DiMethyl-Histone H3-K36 antibody at dilution of 1:200 (40x lens).



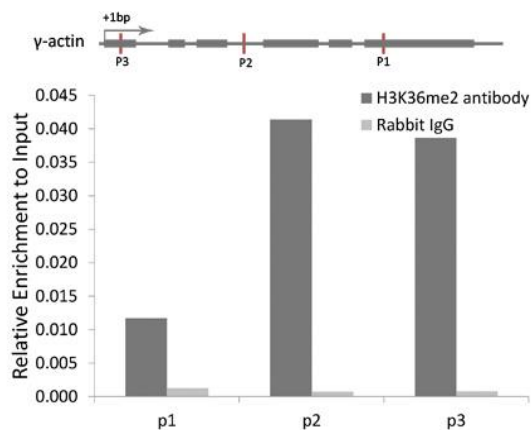
Immunohistochemistry of paraffin-embedded human breast using DiMethyl-Histone H3-K36 antibody at dilution of 1:200 (40x lens).



Immunofluorescence analysis of 293T cells using DiMethyl-Histone H3-K36 antibody. Blue: DAPI for nuclear staining.



Immunohistochemistry of paraffin-embedded rat testis using DiMethyl-Histone H3-K36 antibody at dilution of 1:200 (40x lens).



Chromatin immunoprecipitation analysis of γ -actin gene from 293 cell line, using DiMethyl-Histone H3-K36 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.