
Histone H3K36me3 (H3K36 Trimethyl) Polyclonal Antibody

(Catalog # A-4042)

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 H3K36me3 (H3K36 Trimethyl) 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

Purification

Affinity Purified

Immunogen

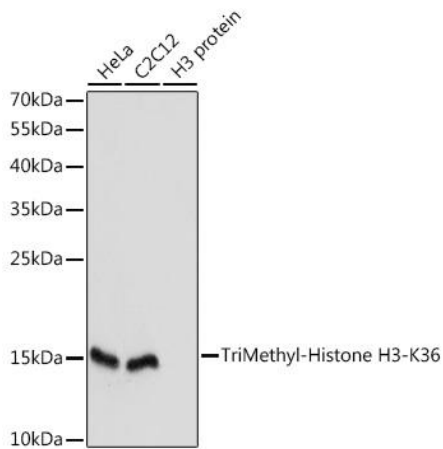
A synthetic peptide corresponding to a sequence within amino acids 1-100 of human histone H3 (NP_003520.1).

Storage

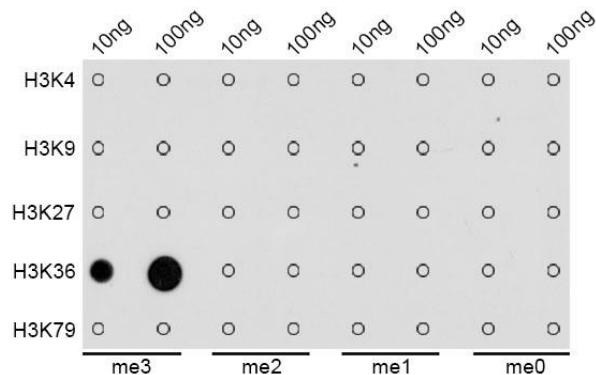
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Application

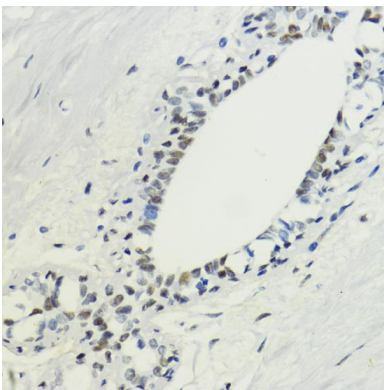
WB, IHC, IF, IP, ChIP, ChIPseq; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:50 - 1:200, IP 1:50 - 1:200, ChIP 1:20 - 1:100, CHIPseq 1:20 - 1:100



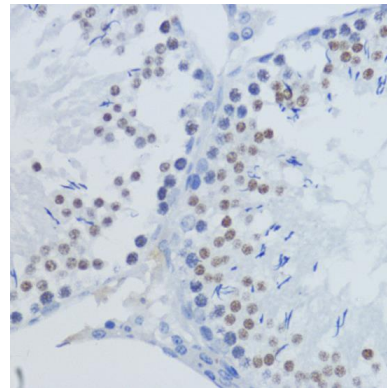
Western blot analysis of extracts of various cell lines, using TriMethyl-Histone H3-K36 antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST.



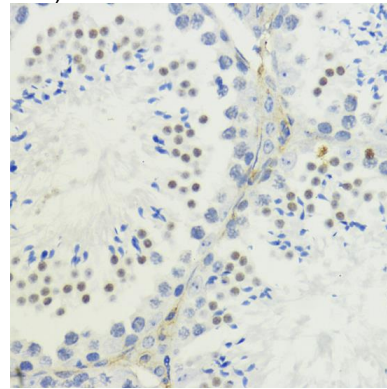
Dot-blot analysis of all sorts of methylation peptides using TriMethyl-Histone H3-K36 antibody.



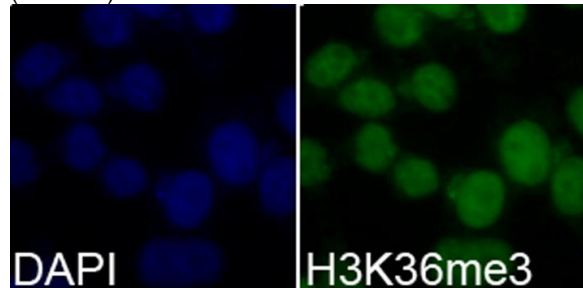
Immunohistochemistry of paraffin-embedded human breast using TriMethyl-Histone H3-K36 antibody at dilution of 1:200 (40x lens).



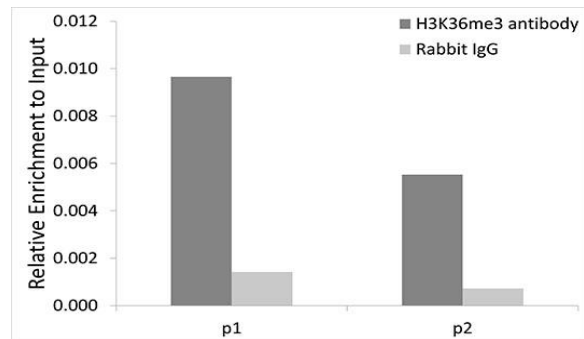
Immunohistochemistry of paraffin-embedded rat testis using TriMethyl-Histone H3-K36 antibody at dilution of 1:200 (40x lens).



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



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



Chromatin immunoprecipitation analysis of extracts of 293T cells, using TriMethyl-Histone H3-K36 antibody and rabbit IgG. P1 and P2 were located on GAPDH gene. The amount of immunoprecipitated DNA was checked by quantitative PCR. Histogram was constructed by the ratios of the immunoprecipitated DNA to the input.