

Histone H3 Methylation Antibody Panel Pack II

- Active Genes

Base Catalog # C10002

PACK CONTENTS

Component	Size	Shipping Temperature	Storage Upon Receipt	Storage Checklist
3K4T Histone H3K4me3 (H3K4 Trimethyl) Polyclonal Antibody	25 µl	4°C	-20°C	
3K36T Histone H3K36me3 (H3K36 Trimethyl) Polyclonal Antibody	25 µl	4°C	-20°C	
3K79T Histone H3K79me3 (H3K79 Trimethyl) Polyclonal Antibody	25 µl	4°C	-20°C	
HGR2 HRP-Goat Anti-Rabbit Secondary Antibody	50 µg	4°C	-20°C	

SHIPPING & STORAGE

This product is shipped on frozen ice packs at 4°C. Upon receipt: (1) Store **3K4T**, **3K36T**, **3K79T**, and **HGR2** at -20°C away from light.

All components of the product are stable for 6 months from the date of shipment, when stored properly.

Histone H3K4me3 (H3K4 Trimethyl) Polyclonal Antibody

Component Cat. #C10002-3K4T

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 H3K4me3 (H3K4 Trimethyl) 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 trimethylated peptide around K4 of human histone H3 (NP_003520.1)

Storage

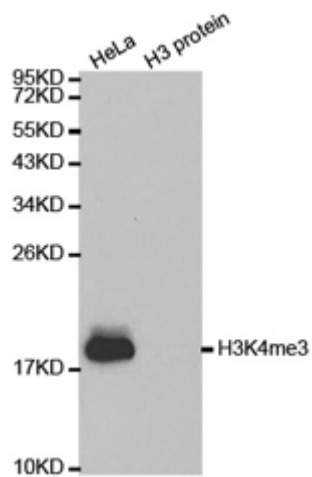
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

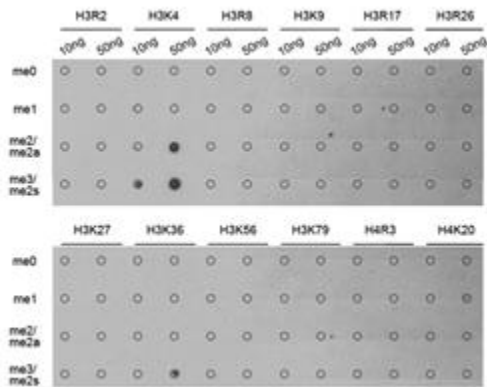
H3K4me3 antibody, H3K4m3 antibody

Application

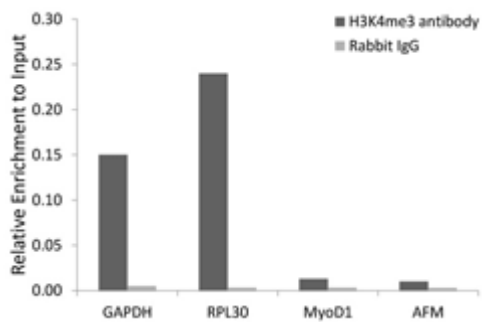
DB, WB, ChIP, ChIP-seq; Recommended dilution, DB 1:500 - 1:2000, WB 1:500 - 1:2000, ChIP 1:20 - 1:100, ChIP-seq 1:20 - 1:100



Western blot analysis of extracts of HeLa cell line and H3 protein expressed in E.coli., using Histone H3K4me3 Polyclonal Antibody.



Dot-blot analysis of all sorts of methylation peptides using H3K4me3 Polyclonal Antibody.



Chromatin immunoprecipitation analysis extracts of 293 cell line, using Histone H3K4me3 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.

Histone H3K36me3 (H3K36 Trimethyl) Polyclonal Antibody

Component Cat. #C10002-3K36T

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

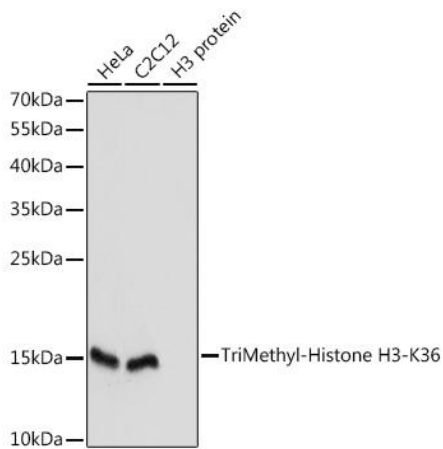
A synthetic peptide corresponding to a sequence within amino acids 1-100 of human histone H3 (NP_003520.1).

Storage

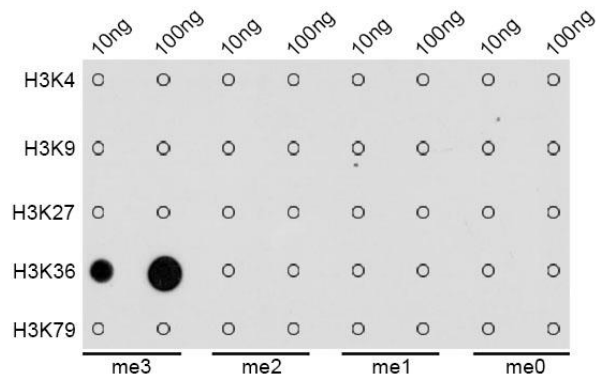
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Application

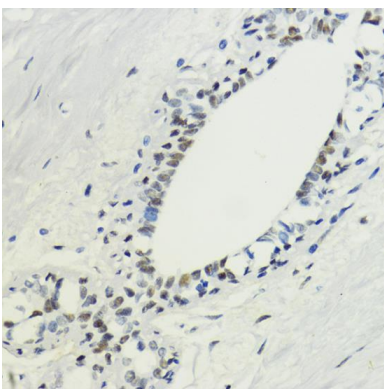
WB, IHC, IF, IP, ChIP, ChIPseq; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:50 - 1:200, IP 1:50 - 1:200, ChIP 1:20 - 1:100, CHIPseq 1:20 - 1:100



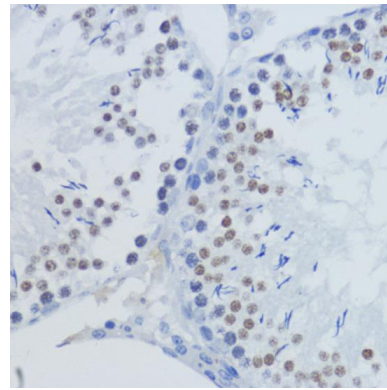
Western blot analysis of extracts of various cell lines, using TriMethyl-Histone H3-K36 antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST.



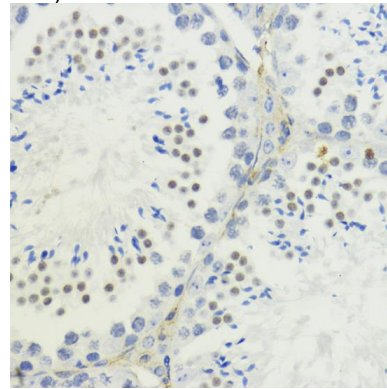
Dot-blot analysis of all sorts of methylation peptides using TriMethyl-Histone H3-K36 antibody.



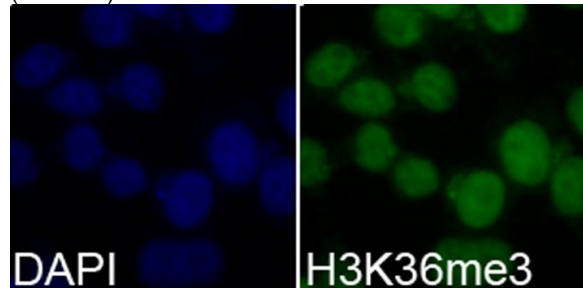
Immunohistochemistry of paraffin-embedded human breast using TriMethyl-Histone H3-K36 antibody at dilution of 1:200 (40x lens).



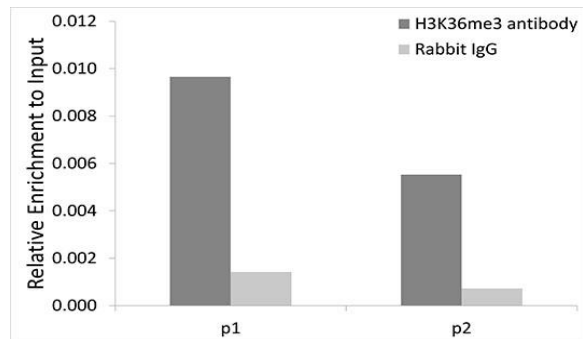
Immunohistochemistry of paraffin-embedded rat testis using TriMethyl-Histone H3-K36 antibody at dilution of 1:200 (40x lens).



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.

Histone H3K79me3 (H3K79 Trimethyl) Polyclonal Antibody

Component Cat. #C10002-3K79T

Description

Histone H3K79me3 (H3K79 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 trimethylated peptide around K79 of human histone H3 (NP_003520.1)

Storage

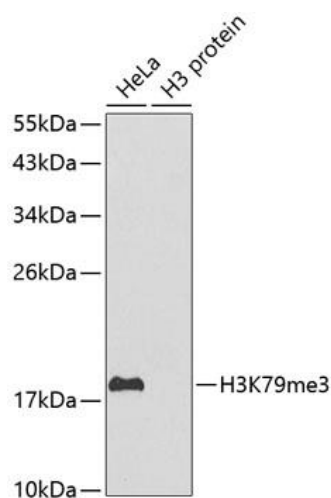
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze.

Alternative Names

HIST3H3, H3/g

Application

DB, WB, IHC, IF, ChIP; Recommended dilution: DB 1:500 - 1:2000, WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:50 - 1:200, ChIP 1:50 - 1:200

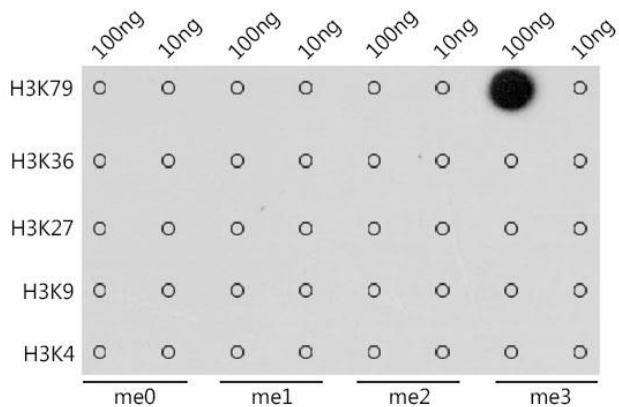


Western blot analysis of extracts of various cell lines, using TriMethyl-Histone H3-K79 antibody.

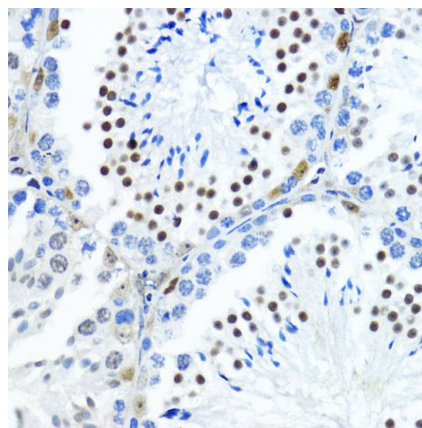
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

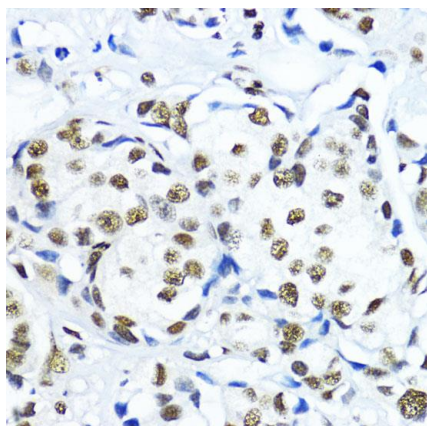
Blocking buffer: 3% nonfat dry milk in TBST.



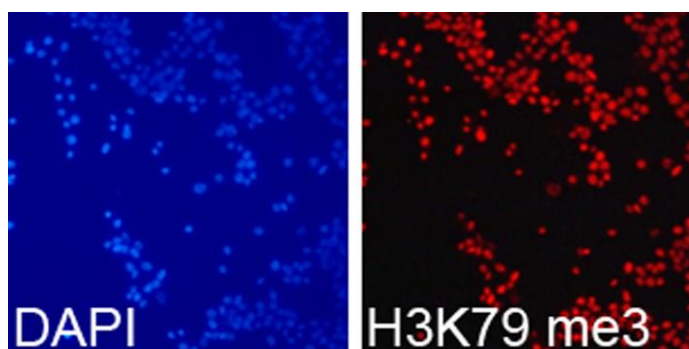
Dot-blot analysis of all sorts of methylation peptides using TriMethyl-Histone H3-K79 antibody at 1:1000 dilution.



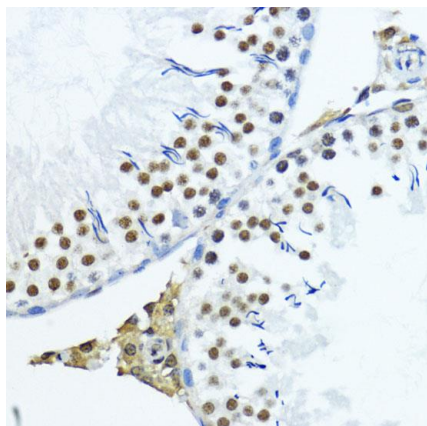
Immunohistochemistry of paraffin-embedded mouse testis using TriMethyl-Histone H3-K79 antibody at dilution of 1:200 (40x lens).



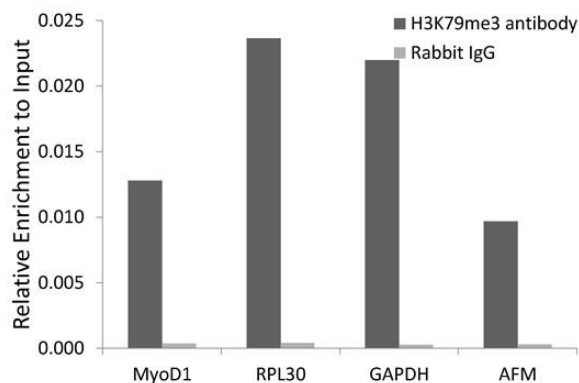
Immunohistochemistry of paraffin-embedded human mammary cancer using TriMethyl-Histone H3-K79 antibody at dilution of 1:200 (40x lens).



Immunofluorescence analysis of 293T cells using TriMethyl-Histone H3-K79 antibody. Blue: DAPI for nuclear staining.



Immunohistochemistry of paraffin-embedded rat testis using TriMethyl-Histone H3-K79 antibody at dilution of 1:200 (40x lens).



Chromatin immunoprecipitation analysis of extracts of 293 cell line, using H3K79me3 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.

HRP- Goat Anti-Rabbit Secondary Antibody

Component Cat. #C10002-HGR2

Description

Goat anti-rabbit IgG recognizes rabbit IgG whole molecule. This secondary antibody was purified using antigen affinity chromatography. The antibody is conjugated with peroxidase.

Antibody Type

Polyclonal Antibody

Purification

Liquid; this product was prepared from monospecific antiserum by immunoaffinity chromatography, followed by solid phase adsorption(s) to remove any unwanted reactivities.

Immunogen

Rabbit IgG whole molecule

Isotype

IgG

Size

350 µg

Formulation

In 10 mM sodium phosphate, 75 mM NaCl, 50% (v/v) glycerol, pH 7.2 , 10 mg/mL BSA (Immunoglobulin and Protease free)

Specificity

Rabbit

Storage

Store at -20°C. Aliquot to avoid repeated freezing and thawing.

Handling Recommendations

The optimal working dilution should be determined by the end user. For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions

Western Blot: 1:1000-1: 10000; Immunohistochemistry:1:100-1:500; Immunofluorescence: 1:100-1:500;
ELISA: 1:2000-1:20000

RELATED PRODUCTS

Histone Modification Antibodies

A-4033	Histone H3K4me3 (H3K4 Trimethyl) Polyclonal Antibody
A-4042	Histone H3K36me3 (H3K36 Trimethyl) Polyclonal Antibody
A-4045	Histone H3K79me3 (H3K79 Trimethyl) Polyclonal Antibody
A12004	HRP-Goat Anti-Rabbit Secondary Antibody

Histone Modification Panel Packs

C10000	Histone H3 Methylation Antibody Panel Pack I – Active Genes
C10001	Histone H3 Methylation Antibody Panel Pack I – Repression Genes
C10002	Histone H3 Methylation Antibody Panel Pack II – Active Genes
C10003	Histone H3 Methylation Antibody Panel Pack II – Repression Genes
C10004	Histone H3 Methylation Antibody Panel Pack III – Active Genes
C10005	Histone H3K4 Methylation Antibody Panel Pack
C10006	Histone H3K9 Methylation Antibody Panel Pack
C10007	Histone H3K27 Methylation Antibody Panel Pack
C10008	Histone H3K36 Methylation Antibody Panel Pack
C10009	Histone H3K79 Methylation Antibody Panel Pack
C10010	Histone H3 Acetylation Antibody Panel Pack I
C10011	Histone H3 Acetylation Antibody Panel Pack II
C10012	Histone H4K20 Methylation Antibody Panel Pack
C10013	Histone H4 Acetylation Antibody Panel Pack
C10014	Histone H3 Phosphorylation Antibody Panel Pack
C10015	Histone H3R2 Methylation Antibody Panel Pack
C10016	Histone H3R8 Methylation Antibody Panel Pack
C10017	Histone H3R17 Methylation Antibody Panel Pack
C10018	Histone H3R26 Methylation Antibody Panel Pack
C10019	Histone H4R3 Methylation Antibody Panel Pack