

## Histone H3K79me3 (H3K79 Trimethyl) Polyclonal Antibody

(Catalog # A-4045)

### 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

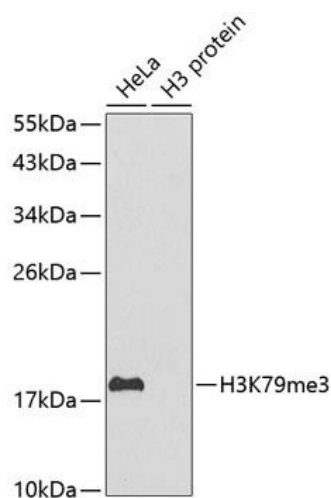
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze.

### 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

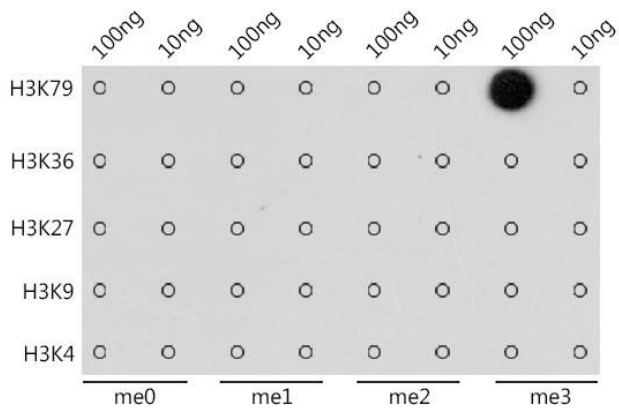


Western blot analysis of extracts of various cell lines, using TriMethyl-Histone H3-K79 antibody.

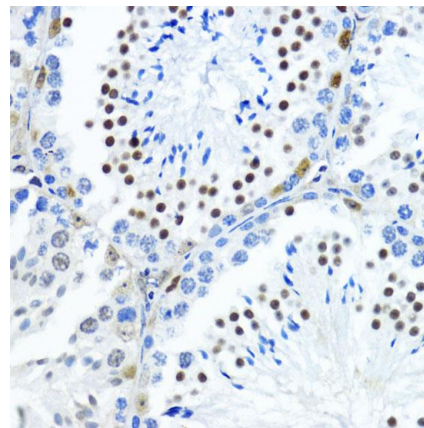
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

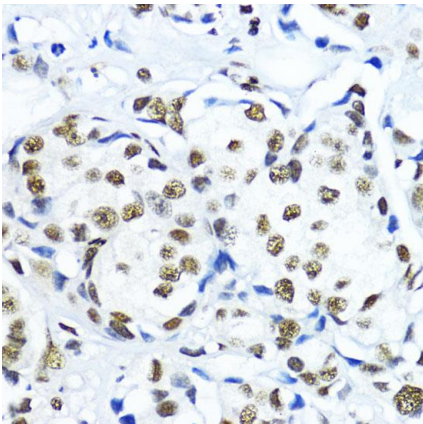
Blocking buffer: 3% nonfat dry milk in TBST.



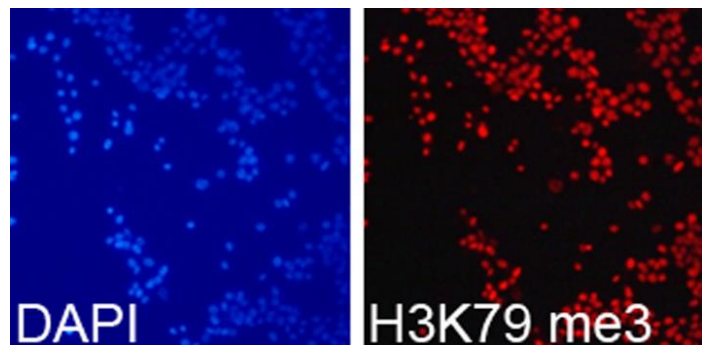
Dot-blot analysis of all sorts of methylation peptides using TriMethyl-Histone H3-K79 antibody at 1:1000 dilution.



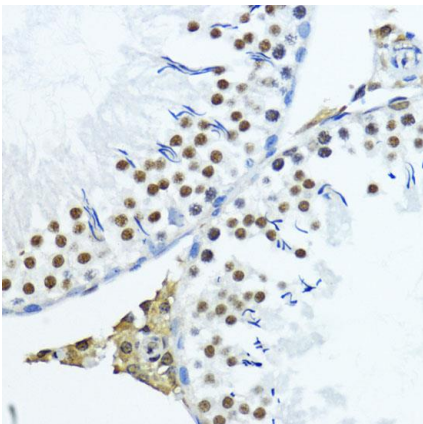
Immunohistochemistry of paraffin-embedded mouse testis using TriMethyl-Histone H3-K79 antibody at dilution of 1:200 (40x lens).



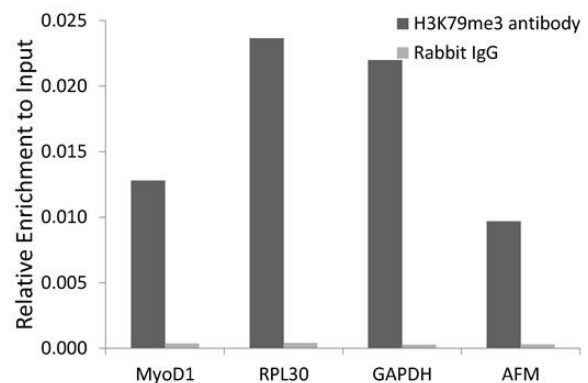
Immunohistochemistry of paraffin-embedded human mammary cancer using TriMethyl-Histone H3-K79 antibody at dilution of 1:200 (40x lens).



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.