

---

## Histone H3R2 Dimethyl Asymmetric (H3R2me2a) Polyclonal Antibody

(Catalog # A-3714)

---

### 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 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 H3R2 Dimethyl Asymmetric (H3R2me2a) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

### Specificity

Human, Mouse, Rat, Broad Range

### Isotype

IgG

### Uniprot ID

Q16695/P68431

### Purification

Affinity Purified

### Immunogen

A synthetic asymmetric dimethylated peptide around R2 of human histone H3 (NP\_003520.1)

### Storage

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

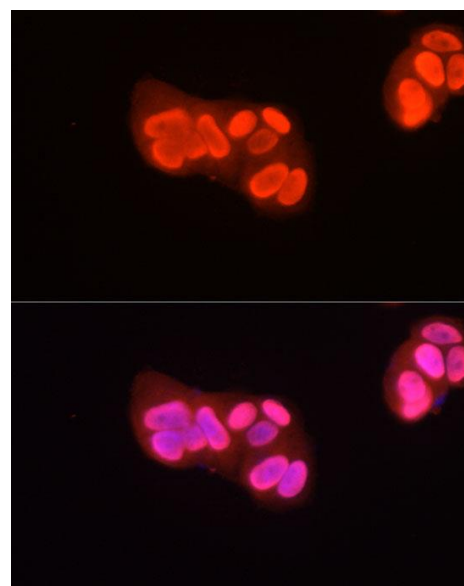
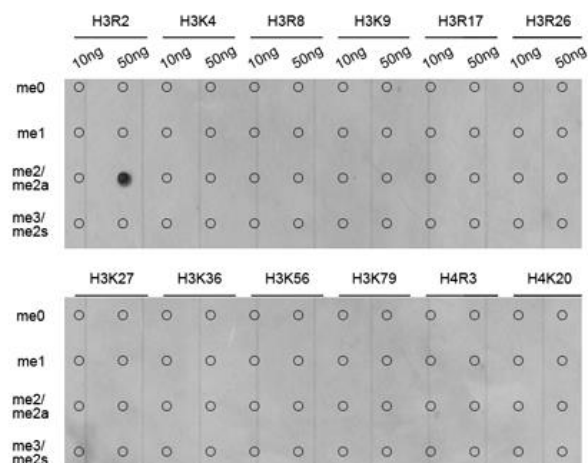
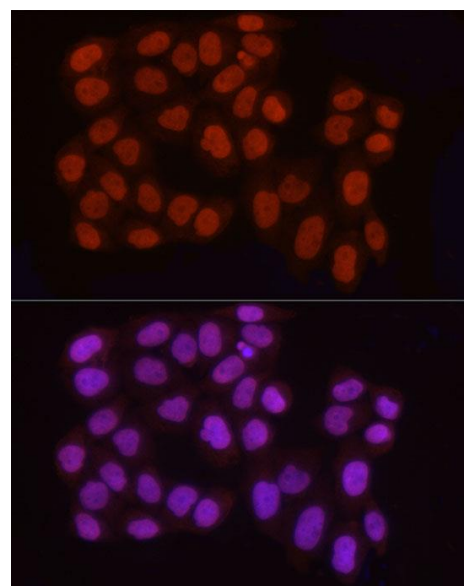
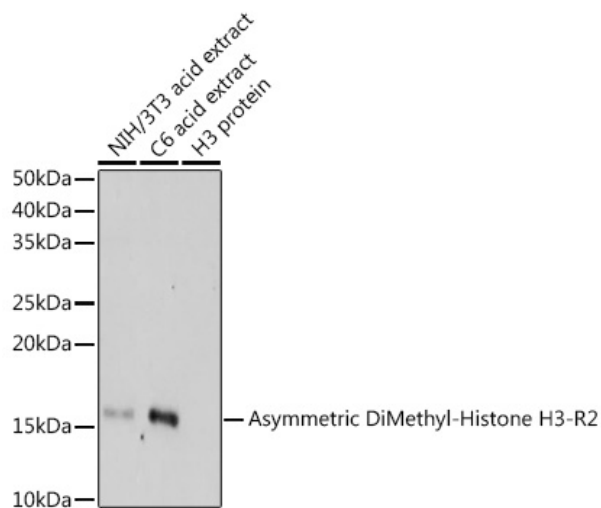
### Alternative Names

H3R2me2a, HIST1H3J, H3/j , H3FJ , Histone H3.1, Histone H3/a, Histone H3/b, Histone H3/c, Histone H3/d, Histone H3/f, Histone H3/h , Histone H3/l, Histone H3/j , Histone H3/k, Histone H3/l, H3 Arginine 2 me2a

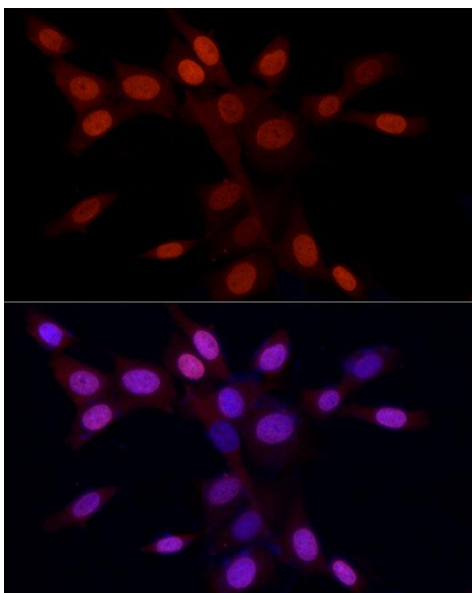
### Application

WB, IF/ICC, ELISA; Recommended dilution: WB 1:100 - 1:500, IF/ICC 1:50 - 1:200, ELISA - recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.

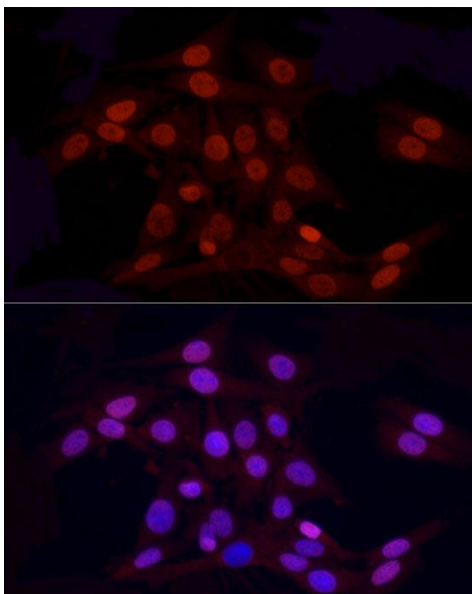
---



Dot-blot analysis of all sorts of methylation peptides using Histone H3R2 Dimethyl Asymmetric (H3R2me2a) Polyclonal Antibody.



Immunofluorescence analysis of NIH/3T3 cells using Histone H3R2 Dimethyl Asymmetric (H3R2me2a) Polyclonal Antibody at dilution of 1:100 (40x lens). Secondary antibody: Cy3-conjugated Goat anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.



Immunofluorescence analysis of PC-12 cells using Histone H3R2 Dimethyl Asymmetric (H3R2me2a) Polyclonal Antibody at dilution of 1:100 (40x lens). Secondary antibody: Cy3-conjugated Goat anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.