

Histone H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody

(Catalog # A-3708)

Description

Histone H4R3 Dimethyl Asymmetric (H4R3me2a) 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

P62805

Purification

Affinity Purified

Immunogen

A synthetic asymmetric dimethylated peptide around R3 of human histone H4 (NP_003539.1)

Storage

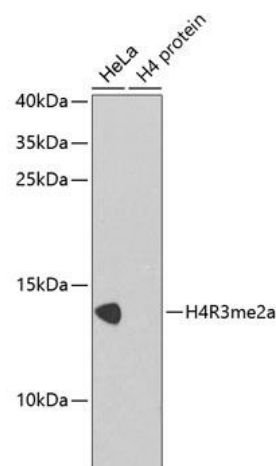
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze.

Alternative Names

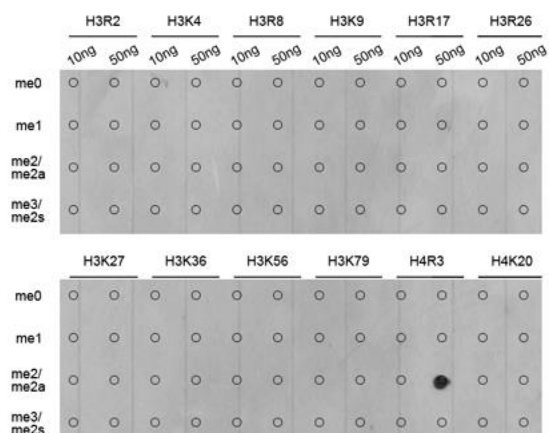
H4R3me2a, H3t, H3.4, H3/g, H3FT, H4 Arginine 3 me2a

Application

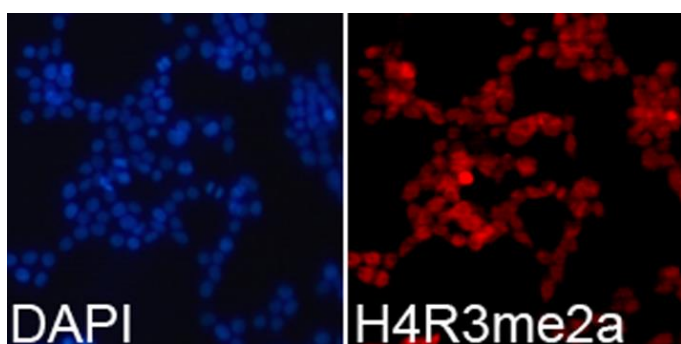
WB, IF, IP; Recommended dilution: WB 1:500 - 1:2000, IF 1:50 - 1:200, IP 1:50 - 1:100



Western blot analysis of extracts of various cell lines, using Asymmetric DiMethyl-Histone H4-R3 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. Lane 1: Mouse liver tissue



Dot-blot analysis of all sorts of methylation peptides using Asymmetric DiMethyl-Histone H4-R3 antibody.



Immunofluorescence analysis of 293T cells using Asymmetric DiMethyl-Histone H4-R3 antibody. Blue: DAPI for nuclear staining.