
Histone H4R3 Dimethyl Symmetric (H4R3me2s) Polyclonal Antibody

(Catalog A-3718)

Background

Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. This structure consists of approximately 146 bp of DNA wrapped around a nucleosome, an 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 H4 family. Transcripts from this gene lack polyA tails; instead, they contain a palindromic termination element. This gene is found in a histone cluster on chromosome 1. This gene is one of four histone genes in the cluster that are duplicated; this record represents the centromeric copy.

Description

Histone H4R3 Dimethyl Symmetric (H4R3me2s) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.09% Sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat, Broad Range

Isotype

IgG

Uniprot ID

P62805

Purification

Affinity Purified

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 1-100 of human Histone H4 (NP_003529.1).

Storage

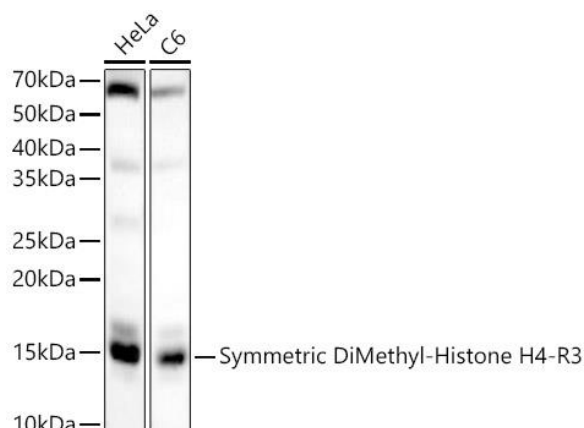
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze.

Alternative Names

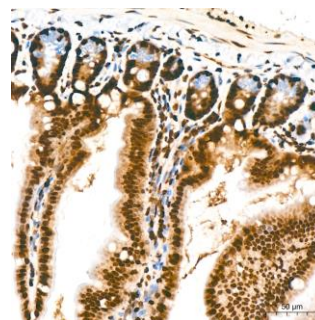
H4R3me2s, H4, H4/n, H4F2, H4FN, FO108, HIST2H4, H4 Arginine 3 me2s

Application

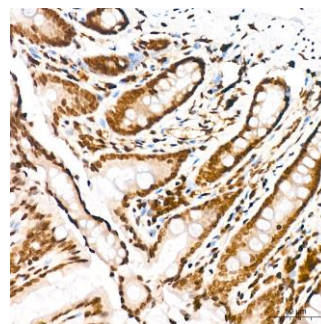
WB, IHC, IF/ICC, DB, ELISA; Recommended dilution: WB 1:500 – 1:2000, IHC 1:50 - 1:200, IF/ICC 1:50 - 1:200, DB 1:500 - 1:2000, ELISA - recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.



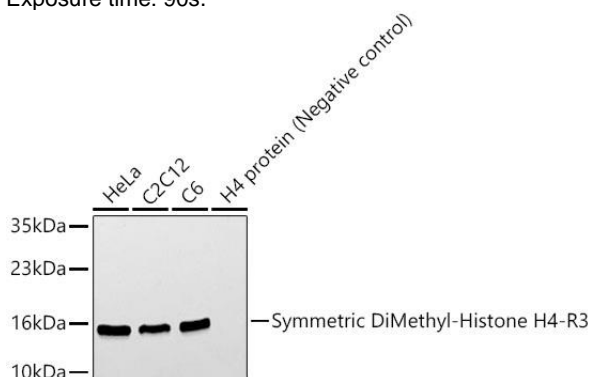
Western blot analysis of various lysates, using H4R3me2s Polyclonal Antibody at 1:2000 dilution.
Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25µg per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 90s.



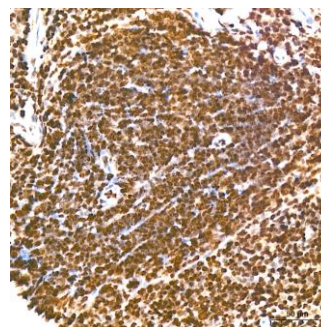
Immunohistochemistry analysis of paraffin-embedded Mouse intestine tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



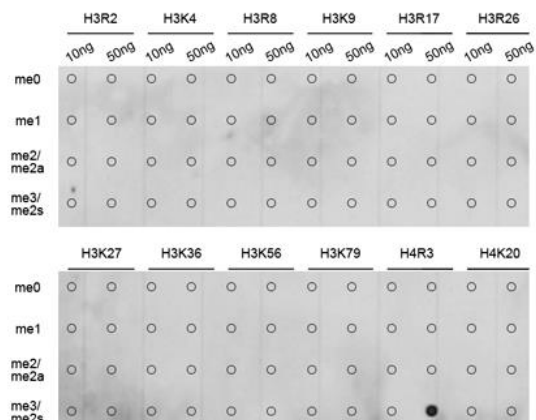
Immunohistochemistry analysis of paraffin-embedded Rat colon tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



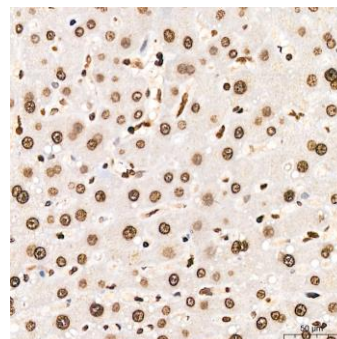
Western blot analysis of various lysates, using H4R3me2s Polyclonal Antibody at 1:400 dilution.
Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25µg per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 60s.



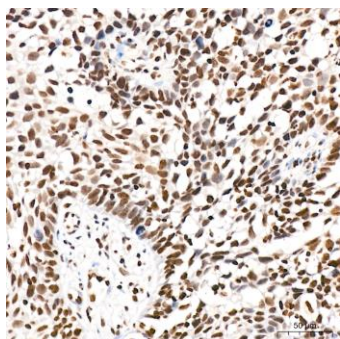
Immunohistochemistry analysis of paraffin-embedded Mouse spleen tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



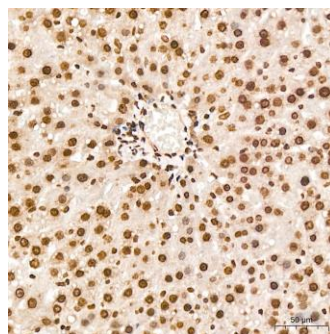
Dot-blot analysis of all sorts of methylation peptides using H4R3me2s Polyclonal Antibody.



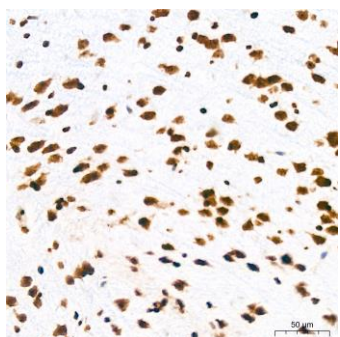
Immunohistochemistry analysis of paraffin-embedded Human liver tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



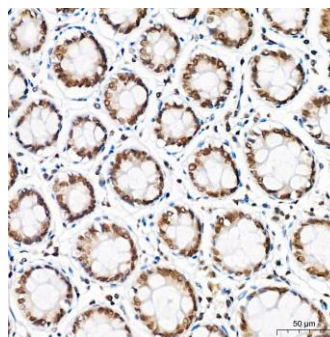
Immunohistochemistry analysis of paraffin-embedded Human cervix cancer tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



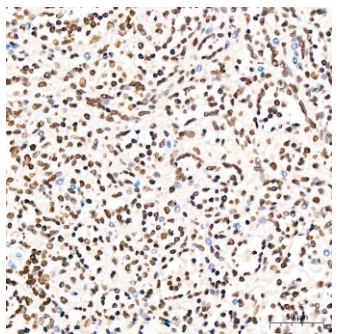
Immunohistochemistry analysis of paraffin-embedded Rat liver tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



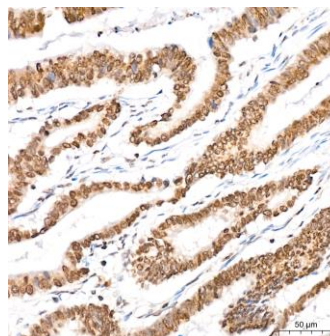
Immunohistochemistry analysis of paraffin-embedded Mouse brain tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



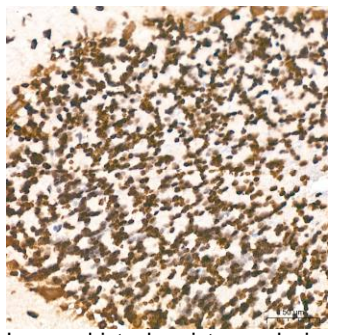
Immunohistochemistry analysis of paraffin-embedded Human colon tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



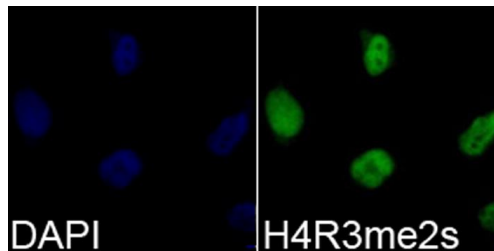
Immunohistochemistry analysis of paraffin-embedded Human spleen tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



Immunohistochemistry analysis of paraffin-embedded Human colon carcinoma tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



Immunohistochemistry analysis of paraffin-embedded Rat brain tissue using H4R3me2s Polyclonal Antibody at a dilution of 1:100 (40x lens). High pressure antigen retrieval was performed with 0.01 M citrate buffer (pH 6.0) prior to IHC staining.



Immunofluorescence analysis of 293T cells using H4R3me2s Polyclonal Antibody. Blue: DAPI for nuclear staining.