
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 0.05% proclin300, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat, Broad Range

Isotype

IgG

Uniprot ID

P62805

Purification

Affinity Purified

Immunogen

A synthetic symmetric dimethylated peptide around R3 of human Histone H4 (NP_003529.1)

Storage

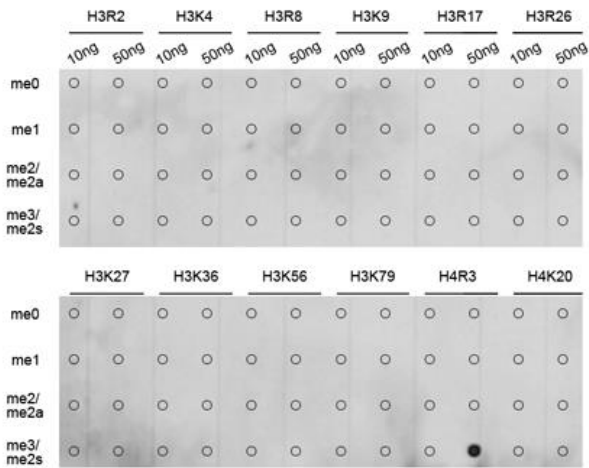
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze.

Alternative Names

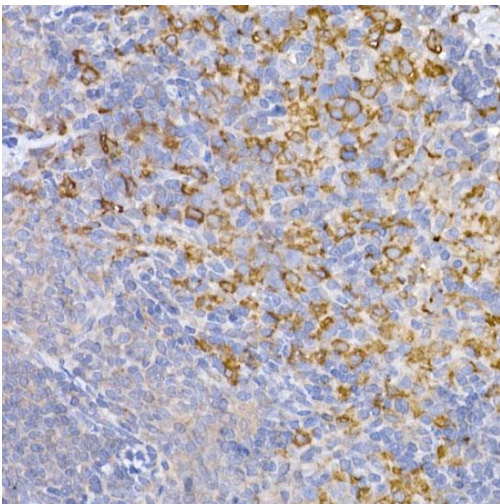
H4R3me2s, H4, H4/n, H4F2, H4FN, FO108, HIST2H4, H4 Arginine 3 me2s

Application

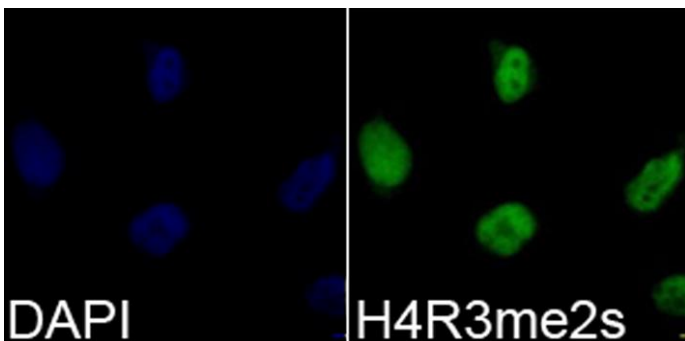
DB, IHC, IF; Recommended dilution: DB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:50 - 1:200



Dot-blot analysis of all sorts of methylation peptides using Histone H4R3 Dimethyl Symmetric (H4R3me2s) Polyclonal Antibody.



Immunohistochemistry of paraffin-embedded mouse brain using Histone H4R3 Dimethyl Symmetric (H4R3me2s) Polyclonal Antibody at dilution of 1:100 (40x lens). Performed high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunofluorescence analysis of 293T cells using Histone H4R3 Dimethyl Symmetric (H4R3me2s) Polyclonal Antibody. Blue: DAPI for nuclear staining.