

Histone H4R3 Methylation Antibody Panel Pack

Base Catalog # C10019

PACK CONTENTS

Component	Size	Shipping Temperature	Storage Upon Receipt	Storage Checklist
4R3M Histone H4R3me1 (H4R3 Monomethyl) Polyclonal Antibody	25 µl	4°C	-20°C	
4R3DA Histone H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody	25 µl	4°C	-20°C	
4R3DS Histone H4R3 Dimethyl Symmetric (H4R3me2s) Polyclonal Antibody	25 µl	4°C	-20°C	
HGR2 HRP-Goat Anti-Rabbit Secondary Antibody	50 µg	4°C	-20°C	

SHIPPING & STORAGE

This product is shipped on frozen ice packs at 4°C. Upon receipt: (1) Store all components at -20°C away from light.

All components are stable for 6 months from the date of shipment, when stored properly.

Histone H4R3 Monomethyl (H4R3me1) Polyclonal Antibody

Component Cat. #C10019-1-4R3M

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 Monomethyl (H4R3me1) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

PBS, pH 7.4, containing 0.02% sodium azide as preservative and 50% glycerol.

Specificity

Human

Isotype

IgG

Uniprot ID

P62805

Purification

Affinity Purified

Immunogen

Synthetic Peptide

Storage

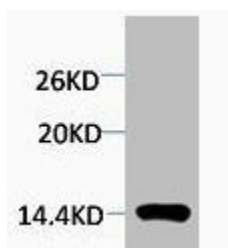
Shipped at 4°C. Aliquot store at -20°C. Avoid repeated freeze / thaw cycles.

Alternative Names

H4R3me1, H4, H4/n, H4F2, H4FN, FO108, HIST2H4, H4 Arginine 3 me1

Application

WB; Recommended dilution: WB: 1:1000-2000



Western blot analysis of extracts from HeLa cells, 1:2000.

Histone H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody

Component Cat. #C10019-1-4R3DA

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 replication-dependent histone that is 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 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

Synthetic Peptide of Human Asymmetric DiMethyl-Histone H4-R3

Storage

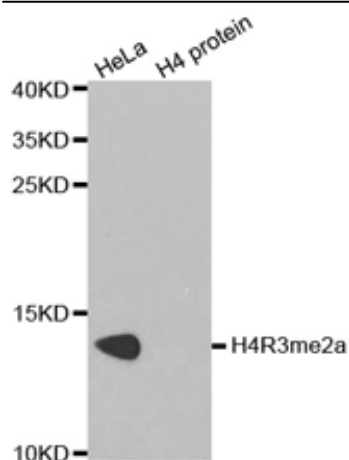
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

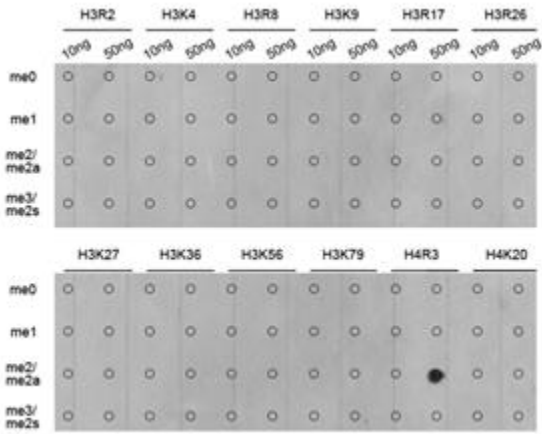
H4R3me2a, H3t, H3.4, H3/g, H3FT, H4 Arginine 3 me2a

Application

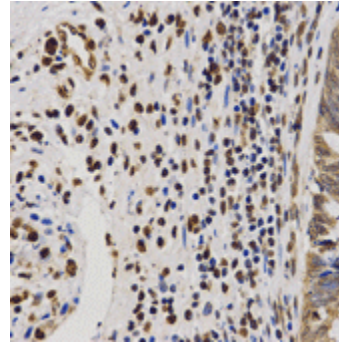
WB, IHC, IF, IP; Recommended dilutions: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:50 - 1:200, IP 1:50 - 1:100



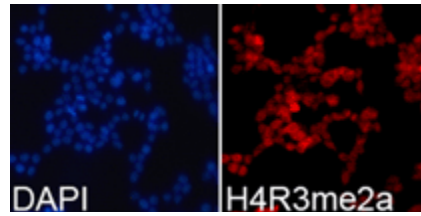
Western blot analysis of Histone H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody.



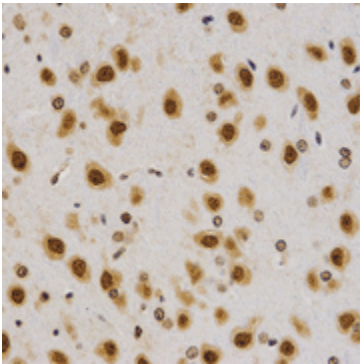
Dot-blot analysis of all sorts of methylation peptides using Histone H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody.



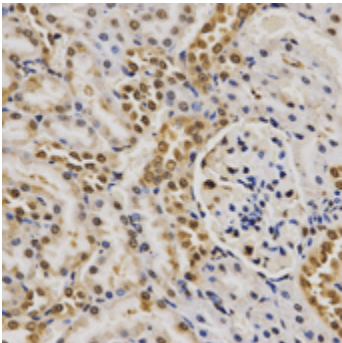
Immunohistochemistry - Histone H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody



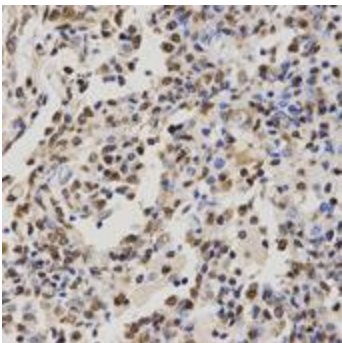
Immunofluorescence analysis -H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody.



Immunohistochemistry - Histone H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody



Immunohistochemistry - Histone H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody



Immunohistochemistry - Histone H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody

Histone H4R3 Dimethyl Symmetric (H4R3me2s) Polyclonal Antibody

Component Cat. #C10019-1-4R3DS

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

Liquid. 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

Synthetic Peptide of Human Symmetric DiMethyl-Histone H4-R3

Storage

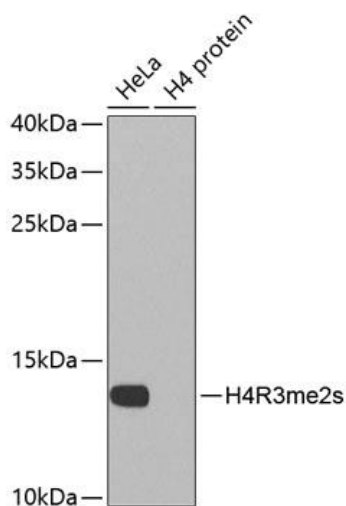
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

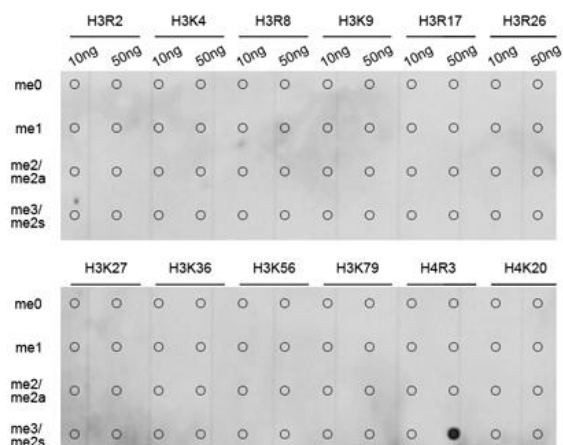
H4R3me2s, H4, H4/n, H4F2, H4FN, FO108, HIST2H4, H4 Arginine 3 me2s

Application

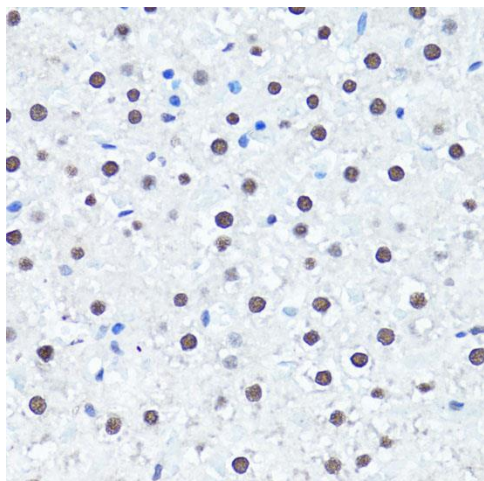
WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:50 - 1:200



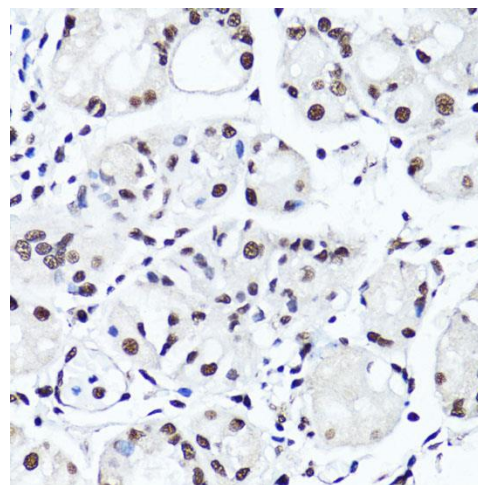
Western blot analysis of extracts of various cell lines, using Symmetric DiMethyl-Histone H4-R3 antibody.
 Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.



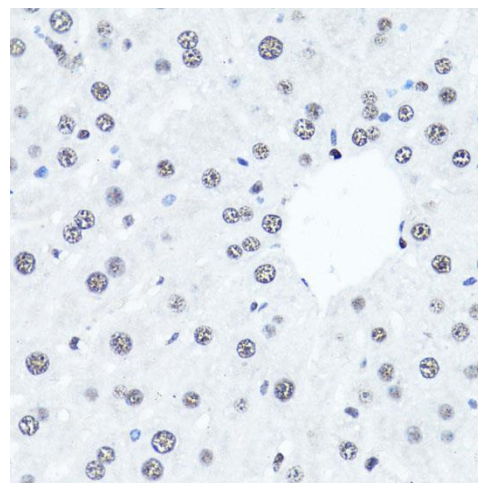
Dot-blot analysis of all sorts of methylation peptides using Symmetric DiMethyl-Histone H4-R3 antibody.



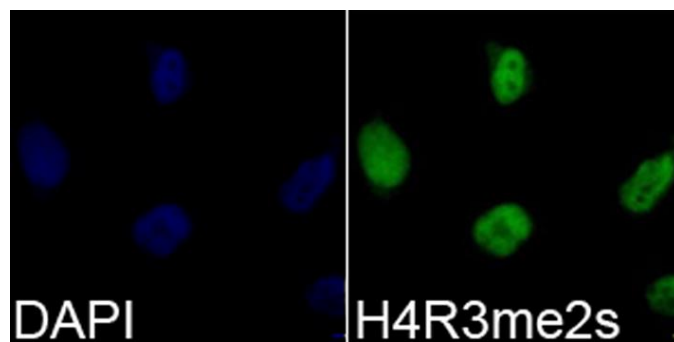
Immunohistochemistry of paraffin-embedded rat liver using Symmetric DiMethyl-Histone H4-R3 antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human stomach using Symmetric DiMethyl-Histone H4-R3 antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded mouse liver using Symmetric DiMethyl-Histone H4-R3 antibody at dilution of 1:100 (40x lens).



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

HRP- Goat Anti-Rabbit Secondary Antibody

Component Cat. #C10019-1-HGR2

Description

Goat anti-rabbit IgG recognizes rabbit IgG whole molecule. This secondary antibody was purified using antigen affinity chromatography. The antibody is conjugated with peroxidase.

Antibody Type

Polyclonal Antibody

Purification

Liquid; this product was prepared from monospecific antiserum by immunoaffinity chromatography, followed by solid phase adsorption(s) to remove any unwanted reactivities.

Immunogen

Rabbit IgG whole molecule

Isotype

IgG

Formulation

In 10 mM sodium phosphate, 75 mM NaCl, 50% (v/v) glycerol, pH 7.2.

Specificity

Rabbit

Storage

Store at -20°C. Aliquot to avoid repeated freezing and thawing.

Handling Recommendations

The optimal working dilution should be determined by the end user. For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions

Western Blot: 1:1000-1: 10000; Immunohistochemistry: 1:100-1:500; Immunofluorescence: 1:100-1:500;
ELISA: 1:2000-1:20000

RELATED PRODUCTS

Histone Modification Antibodies

A-3717	Histone H4R3 Monomethyl (H4R3me1) Polyclonal Antibody
A-3708	Histone H4R3 Dimethyl Asymmetric (H4R3me2a) Polyclonal Antibody
A-3718	Histone H4R3 Dimethyl Symmetric (H4R3me2s) Polyclonal Antibody
A12004	HRP-Goat Anti-Rabbit Secondary Antibody

Histone Modification Panel Packs

C10000	Histone H3 Methylation Antibody Panel Pack I – Active Genes
C10001	Histone H3 Methylation Antibody Panel Pack I – Repression Genes
C10002	Histone H3 Methylation Antibody Panel Pack II – Active Genes
C10003	Histone H3 Methylation Antibody Panel Pack II – Repression Genes
C10004	Histone H3 Methylation Antibody Panel Pack III – Active Genes
C10005	Histone H3K4 Methylation Antibody Panel Pack
C10006	Histone H3K9 Methylation Antibody Panel Pack
C10007	Histone H3K27 Methylation Antibody Panel Pack
C10008	Histone H3K36 Methylation Antibody Panel Pack
C10009	Histone H3K79 Methylation Antibody Panel Pack
C10010	Histone H3 Acetylation Antibody Panel Pack I
C10011	Histone H3 Acetylation Antibody Panel Pack II
C10012	Histone H4K20 Methylation Antibody Panel Pack
C10013	Histone H4 Acetylation Antibody Panel Pack
C10014	Histone H3 Phosphorylation Antibody Panel Pack
C10015	Histone H3R2 Methylation Antibody Panel Pack
C10016	Histone H3R8 Methylation Antibody Panel Pack
C10017	Histone H3R17 Methylation Antibody Panel Pack
C10018	Histone H3R26 Methylation Antibody Panel Pack
C10019	Histone H4R3 Methylation Antibody Panel Pack