
Histone H4K20me1 (H4K20 Monomethyl) Polyclonal Antibody

(Catalog # A-4046)

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 H4K20me1 (H4K20 Monomethyl) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Broad Range, Mouse, Rat, Human

Isotype

IgG

Uniprot ID

P62805

Purification

Affinity Purified

Immunogen

A synthetic monomethylated peptide around K20 of human Histone H4 (NP_003529.1).

Storage

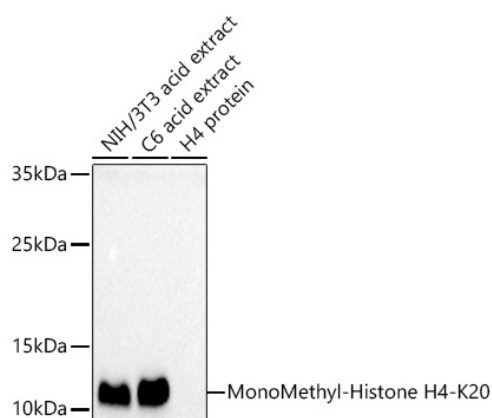
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

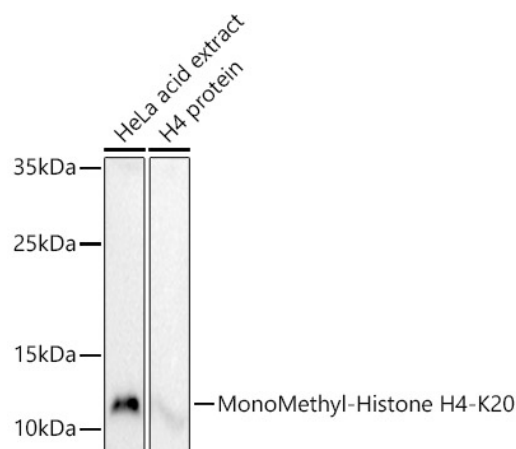
H4; H4/n; H4F2; H4FN; FO108; HIST2H4; H4K20me1 antibody; H4K20m1 antibody

Applications

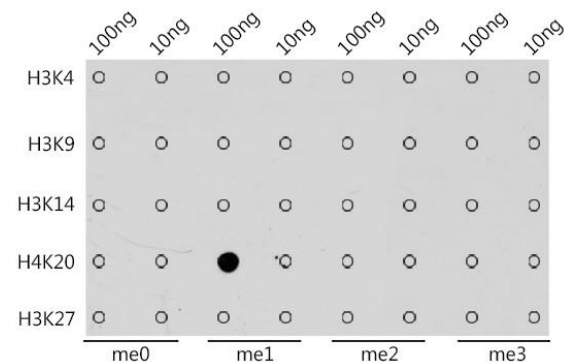
WB, DB, IHC, IF/ICC, ELISA; Recommended dilution: WB 1:500 - 1:1000, DB 1:500 - 1:2000, IHC 1:50 - 1:200, IF/ICC 1:50 - 1:200, ELISA - recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.



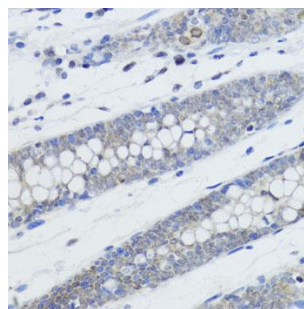
Western blot analysis of various lysates using Histone H4K20me1 (H4K20 Monomethyl) Polyclonal Antibody at 1:1000 dilution.
 Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25µg per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.
 Exposure time: 10s.



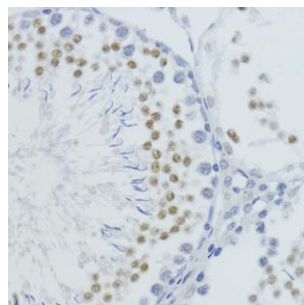
Western blot analysis of various lysates using Histone H4K20me1 (H4K20 Monomethyl) Polyclonal Antibody at 1:1000 dilution.
 Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25µg per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.
 Exposure time: 90s.



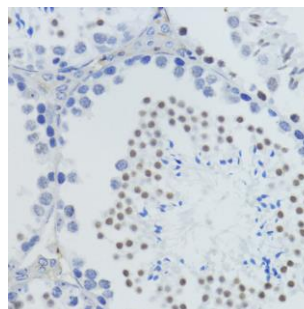
Dot-blot analysis of all sorts of methylation peptides using Histone H4K20me1 (H4K20 Monomethyl) Polyclonal Antibody.



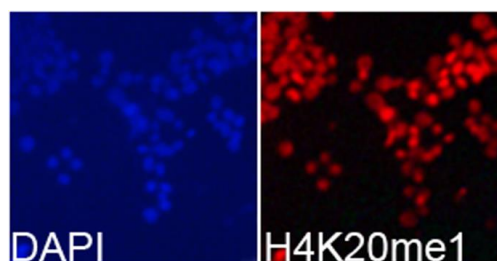
Immunohistochemistry analysis of paraffin-embedded Human colon using Histone H4K20me1 (H4K20 Monomethyl) Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



Immunohistochemistry analysis of paraffin-embedded Rat testis using Histone H4K20me1 (H4K20 Monomethyl) Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



Immunohistochemistry analysis of paraffin-embedded Mouse testis using Histone H4K20me1 (H4K20 Monomethyl) Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



Immunofluorescence analysis of 293T cells using Histone H4K20me1 (H4K20 Monomethyl) Polyclonal Antibody. Blue: DAPI for nuclear staining.