
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

PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Broad Range, Mouse, Rat, Human

Isotype

IgG

Uniprot ID

P62805

Purification

Affinity Purified

Immunogen

Synthetic Peptide of Human MonoMethyl-Histone H4-K20

Storage

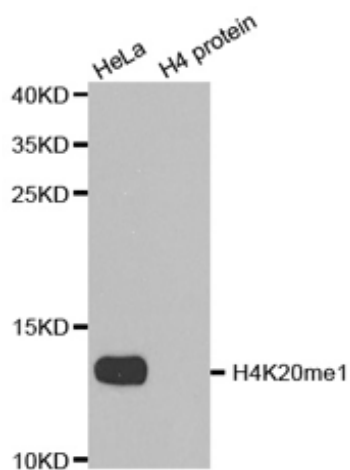
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

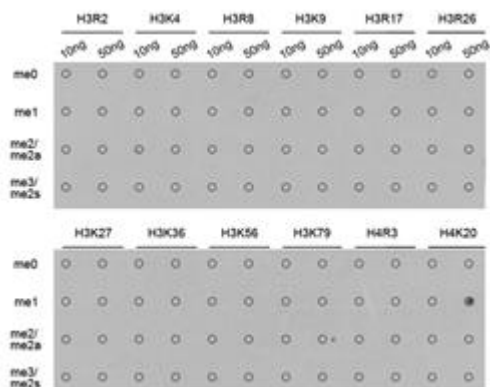
H4; H4/n; H4F2; H4FN; FO108; HIST2H4; H4K20me1 antibody; H4K20m1 antibody

Application

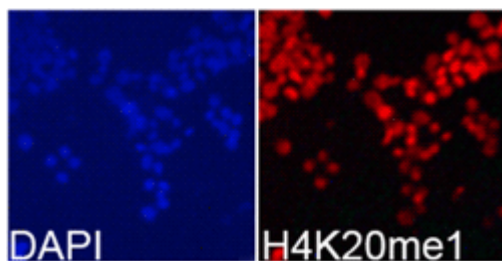
WB, IHC, IF, IP, ChIP, ChIPseq; Recommended dilution: WB 1:500 - 1:2000 IHC 1:50 - 1:200 IF 1:50 - 1:200 IP 1:50 - 1:200 ChIP 1:20 - 1:100 CHIPseq 1:20 - 1:100



Western blot analysis of extracts of HeLa cell line and H4 protein expressed in E.coli., using H4K20me1 Polyclonal Antibody.



Dot-blot analysis of all sorts of methylation peptides using H4K20me1 Polyclonal Antibody.



Immunofluorescence analysis of 293T cell using H4K20me1 Polyclonal Antibody. Blue: DAPI for nuclear staining.