
Histone H3.1K9me2 (H3.1K9 Dimethyl) Monoclonal Antibody [RMC418M]

(Catalog # A68435)

Background

Core component of nucleosome. Nucleosomes wrap and compact DNA into chromatin, limiting DNA accessibility to the cellular machineries which require DNA as a template. Histones thereby play a central role in transcription regulation, DNA repair, DNA replication and chromosomal stability. DNA accessibility is regulated via a complex set of post-translational modifications of histones, also called histone code, and nucleosome remodeling.

Description

Histone H3.1K9me2 (H3.1K9 Dimethyl) Monoclonal Antibody, clone RMC418M. Unconjugated. Raised in: Rabbit.

Concentration

Inquire

Formulation

Liquid. Rabbit IgG in phosphate buffered saline, pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Specificity

Human

Isotype

IgG

Uniprot ID

P68431

Purification

Affinity Purified

Immunogen

A synthesized peptide

Storage

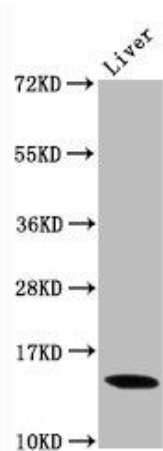
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Histone H3.1, Histone H3/a, Histone H3/b, Histone H3/c, Histone H3/d, Histone H3/f, Histone H3/h, Histone H3/i, Histone H3/j, Histone H3/k, Histone H3/l, HIST1H3A, H3FA, AND, HIST1H3B, H3FL, AND, HIST1H3C, H3FC, AND, HIST1H3D, H3FB, AND, HIST1H3E, H3FD, AND, HIST1H3F, H3FI, AND, HIST1H3G, H3FH, AND, HIST1H3H, H3FK, AND, HIST1H3I, H3FF, AND, HIST1H3J, H3FJ

Application

ELISA, WB, IF; Recommended dilution: WB: 1:500-1:5000, IF: 1:30-1:200



Western Blot

Positive WB detected in: Mouse liver tissue

All lanes: Histone H3.1K9me2 (H3.1K9 Dimethyl) Monoclonal

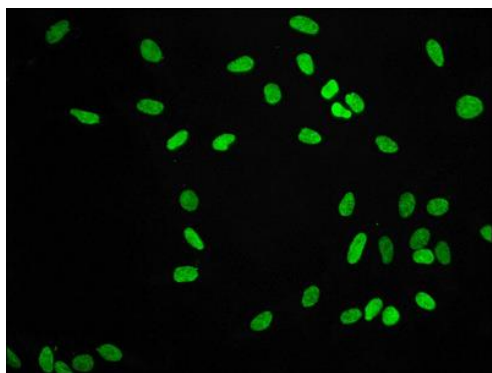
Antibody [RMC418M] at 0.9 ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 15 KDa

Observed band size: 15 KDa



Immunofluorescence staining of HeLa cells with Histone H3.1K9me2 (H3.1K9 Dimethyl) Monoclonal Antibody [RMC418M] at 1:56, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).