

Histone H2AB/E Monoclonal Antibody [RMC385C]

(Catalog # A68424)

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 H2AB/E Monoclonal Antibody, clone RMC385C. 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

P04908

Purification

Affinity Purified

Immunogen

A synthesized peptide

Storage

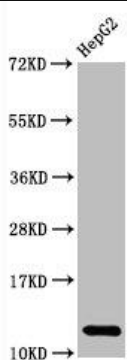
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Histone H2A type 1-B/E, Histone H2A.2, Histone H2A/a, Histone H2A/m, HIST1H2AB, H2AFM, AND, HIST1H2AE, H2AFA

Application

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



Western Blot

Positive WB detected in: HepG2 whole cell lysate

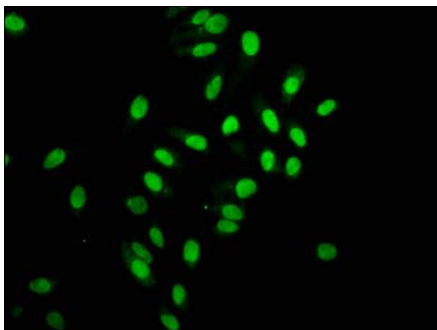
All lanes: Histone H2AB/E Monoclonal Antibody [RMC385C] at 2.7 ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 13 kDa

Observed band size: 13 kDa



Immunofluorescence staining of HeLa cells with Histone H2AB/E Monoclonal Antibody [RMC385C] at 1:168, 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).