

Histone H2A.2K9ac (Acetyl H2A.2K9) Recombinant Monoclonal Antibody [3H5]

(Catalog #A73825)

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 H2A.2K9ac (Acetyl H2A.2K9) Recombinant Monoclonal Antibody [3H5]. Unconjugated. Raised in: HEK293F Cell.

Formulation

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

Specificity

Human

Isotype

Rabbit IgG

Uniprot ID

P04908

Purification

Affinity Chromatography

Immunogen

A synthesized peptide

Storage

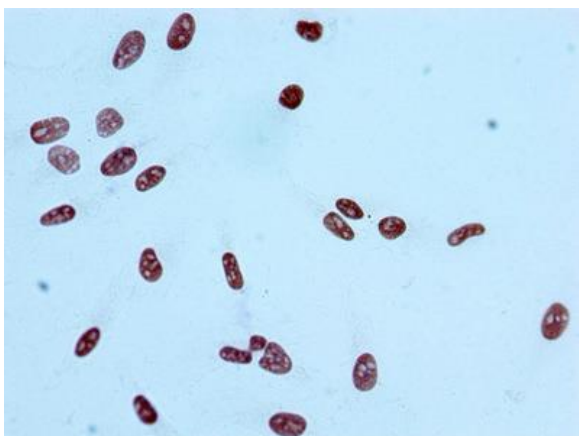
Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

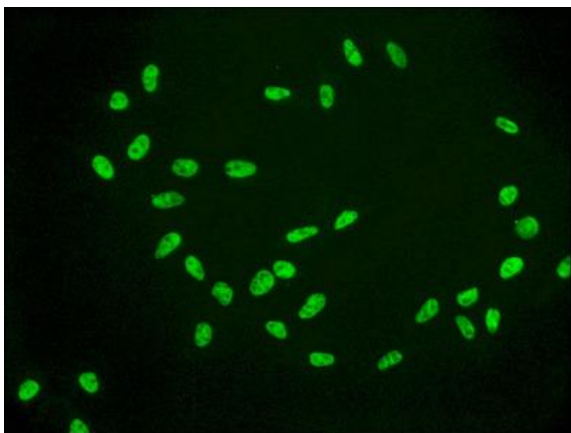
Histone H2A type 1-B/E, Histone H2A.2, Histone H2A/a, Histone H2A/m, HIST1H2AB, H2AFM, HIST1H2AE, H2AFA, Acetyl-Histone H2A type 1-B/E (K9)

Application

ELISA, ICC, IF; Recommended dilution: ICC:1:50-1:500, IF:1:30-1:200



Immunocytochemistry analysis of Histone H2A.2K9ac rMAb diluted at 1:100 and staining in HeLa cells performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Immunofluorescence staining of HeLa cells (treated by 15mM sodium butyrate for 30min) with Histone H2A.2K9ac rMAb 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).