

## Histone H3K9ac (Acetyl H3K9) Monoclonal Antibody [4E9B11]

(Catalog # A-4054)

### Background

Histone H3- along with H2A, H2B and H4- is involved in the structure of chromatin in eukaryotic cells. Histone H3 can undergo several different types of epigenetic modifications that influence cellular processes. These modifications including acetylation, phosphorylation, methylation, ubiquitination, and ADP-ribosylation occur on the N-terminal tail domains of histone H3, which results in remodeling of the nucleosome structure into an open conformation more accessible to transcription complexes. In most species, histone H3 is primarily acetylated at lysine 9, 14, 18, and 23.

### Description

Mouse monoclonal antibody raised against a synthetic peptide corresponding to the N-terminus of histone H3 acetylated on K9

### Concentration

1 mg/ml

### Purification

Protein A purified

### Specificity

Detects histone H3 only when acetylated at K9 in mouse, rat, and human

### Isotype

IgG

### Formulation

PBS (pH 7.5), 0.03% NaN<sub>3</sub>

### Storage

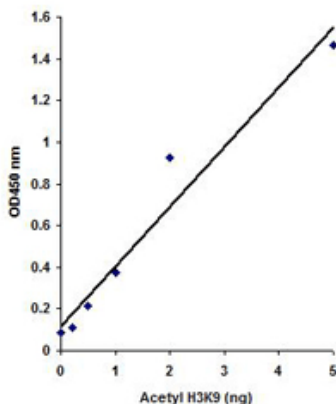
Store at -20°C

### Alternative Names

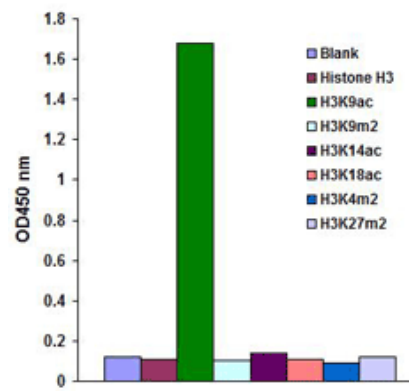
H3K9ac antibody, H3K9a antibody

### Application

WB: 1:200 – 1:1000, IHC: 1:100 – 1:500, ELISA: 1:1000 – 1:2000



High sensitivity of Histone H3K9ac (Acetyl H3K9) Monoclonal Antibody [4E9B11] in detecting histone H3 acetylated at K9.



High specificity of Histone H3K9ac (Acetyl H3K9) Monoclonal Antibody [4E9B11] in detecting histone H3 acetylated at K9.