

AbX™

# Histone H3 Rabbit pAb

Catalog Number A004-100UL



ARBOR  
ASSAYS

## FEATURES

- Reacts with modified Histone H3
- Supplied as a PBS solution
- Applications include Western blotting and ChIP

## INTRODUCTION

Chromosomal surfaces are ornamented with a variety of posttranslational modifications of histones, which are required for the regulation of many of the DNA-templated processes. Histone H3 is one of the five main histone proteins involved in the structure of chromatin in eukaryotic cells. Featuring a main globular domain and a long N-terminus tail, H3 is involved with the 'beads on a string' structure of the nucleosome. Histone H3 is the most extensively post-translationally modified of the five histones. The term "Histone H3" alone is purposely ambiguous in that it does not distinguish between sequence variants or modification state. Histone H3 is an important protein in the emerging field of epigenetics, where its sequence variants and variable modification states are thought to play a role in the regulation of genes. For example, histone methylation either on their arginine and/or lysine residues, yields a more diverse array of cellular functions.

## FORM

AbX™ C-Terminal Histone H3 Rabbit Antibody is produced as diluted rabbit serum.

## IMMUNOGEN

C-Terminal peptide of Histone H3 which should be non-modified.

## CROSS REACTIVITY

Peptide immunogen sequence is highly conserved. Expected to react with Histone H3 from yeast to mammals.

## BUFFER COMPOSITION

Phosphate Buffered Saline at pH 7.2 containing 0.1% Tween 20 and 0.09% Kathon preservative

## STORAGE

Short Term: 4°C. Extended: Aliquot and freeze at -20°C

## USES

Western blotting and Immunoprecipitation

## SUGGESTED DILUTION

Western blotting: Suggested dilution, 1:500-1:1,500  
ChIP: 50 µL

**FOR RESEARCH USE ONLY**

info@ArborAssays.com • www.ArborAssays.com

**WESTERN BLOTTING**

5  $\mu$ L of acid treated HeLa extracts or recombinant Histone H3 were treated with a 1:500 dilution of antibody A004-100UL prior to visualization.

**Related Products**

DetectX™ Histone Demethylase Fluorescent Activity Kit  
Catalog Number K010-F1

AbX™ Antibodies to LSDI and Unmodified, Monomethylated,  
Dimethylated and Trimethylated Lysine<sup>4</sup> Histone H3

Catalog Numbers A003, A005, A006, and A007