

# Histone H3 Polyclonal Antibody

(Catalog # A-1112)

## Background

The primary building block of chromatin is the nucleosome which is made up of DNA wound around eight core histone proteins (two each of H2A, H2B, H3, and H4). Histones thereby play a central role in transcription regulation, DNA repair, DNA replication, and chromosomal stability. DNA accessibility is regulated through an intricate series of post-translational modifications of histones, known as the histone code, and nucleosome remodeling. Histone H3 is mainly acetylated at Lys9, 14, 18, 23, 27, and 56. Acetylation of H3 at Lys9 appears to have a dominant role in histone deposition and chromatin assembly in some organisms.

## Description

Histone H3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

PBS, pH 7.4, containing 0.02% sodium azide as Preservative and 50% Glycerol.

## Immunogen

Recombinant Protein

## Purification

Affinity Purified

## Specificity

Human

## Isotype

IgG

## Uniprot ID

P68431, Q71DI3, P84243

## Alternative Names

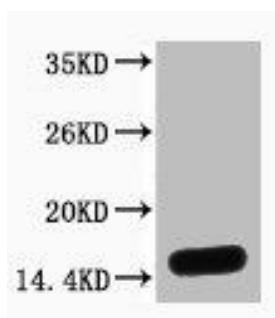
Anti-histone H3, H3 histone antibody

## Storage

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

## Application

WB, ELISA; Recommended dilution: WB: 1:500-1:2000, ELISA 1:40000



Western blot analysis of HeLa, diluted at 1:2000 using the Histone H3 Polyclonal Antibody.