

# Histone H3K14me3 (H3K14 Trimethyl) Polyclonal Antibody

(Catalog # A-3723)

## Background

Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. Nucleosomes consist of approximately 146 bp of DNA wrapped around a histone octamer composed of pairs of each of the four core histones (H2A, H2B, H3, and H4). The chromatin fiber is further compacted through the interaction of a linker histone, H1, with the DNA between the nucleosomes to form higher order chromatin structures. This gene is intronless and encodes a member of the histone H3 family. Transcripts from this gene lack polyA tails; instead, they contain a palindromic termination element. This gene is located separately from the other H3 genes that are in the histone gene cluster on chromosome 6p22-p21. 3.

## Description

Histone H3K14me3 (H3K14 Trimethyl) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

Liquid. PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

## Specificity

Human, Mouse, Rat, Broad Range

## Isotype

IgG

## Uniprot ID

Q16695

## Purification

Affinity Purified

## Immunogen

Synthetic Peptide of Human TriMethyl-Histone H3-K14

## Storage

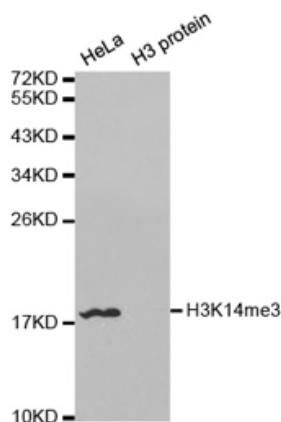
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

## Alternative Names

H3t, H3.4, H3/g, H3FT

## Application

WB, IHC, IF, IP, ChIP, ChIPseq; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:50 - 1:200, IP 1:50 - 1:200, ChIP 1:20 - 1:100, CHIPseq 1:20 - 1:100



Western blot - Histone H3K14 Trimethyl Polyclonal Antibody