
Phospho Histone H3.3 (Ser32) Polyclonal Antibody

(Catalog # A-7700)

Background

Histones are proteins found in eukaryotic cell nuclei that package and order the DNA into structural units called nucleosomes. Nucleosomes consist of about 146-147 bp of DNA wrapped around an octamer of histone proteins (histone 2A, histone 2B, histone 3, and histone 4). The N-terminal tails of histones protrude from the globular nucleosome core and can undergo several different types of epigenetic modifications that influence cellular processes. The interaction of linker histone, H1, with DNA between nucleosomes facilitates the compaction of chromatin into higher-order structures. Unlike most histone genes, this gene contains introns and its mRNA is polyadenylated. The protein encoded is a replication-independent member of the histone H3 family. Histone variant H3.3 is typically enriched in active chromatin.

Description

Phospho Histone H3.3 (Ser32) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P84243

Purification

Affinity Purified

Storage

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

Alternative Names

H3F3, H3.3A; H3.3Ser32; H3.3Ser32ph

Application

WB; Recommended dilution: WB 1:500 - 1:2000
