

H3F3A (Ab-79) Polyclonal Antibody

(Catalog # A70375)

Background

Variant histone H3 which replaces conventional H3 in a wide range of nucleosomes in active genes. Constitutes the predominant form of histone H3 in non-dividing cells and is incorporated into chromatin independently of DNA synthesis. Deposited at sites of nucleosomal displacement throughout transcribed genes, suggesting that it represents an epigenetic imprint of transcriptionally active chromatin. Nucleosomes wrap and compact DNA into chromatin, limiting DNA accessibility to the cellular machineries which require DNA as a template. Histones thereby play a central role in transcription regulation, DNA repair, DNA replication and chromosomal stability. DNA accessibility is regulated via a complex set of post-translational modifications of histones, also called histone code, and nucleosome remodeling.

Description

H3F3A (Ab-79) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300. 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human, Rat

Isotype

IgG

Uniprot ID

P84243

Purification

Affinity purified

Immunogen

Peptide sequence around site of Lys (79) derived from Human Histone H3.3

Storage

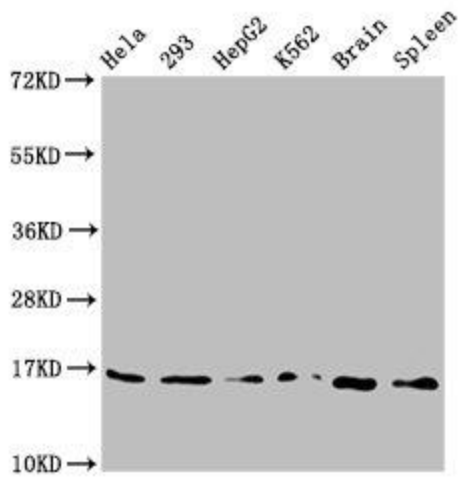
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Histone H3.3, H3F3A, H3.3A, H3F3, PP781, AND, H3F3B, H3.3B

Application

ELISA, WB, ChIP; Recommended dilution: WB:1:200-1:2000



Western Blot

Positive WB detected in: HeLa whole cell lysate, 293 whole cell lysate, HepG2 whole cell lysate, K562 whole cell lysate, Rat brain tissue, Rat spleen tissue

All lanes: H3F3A antibody at 0.73ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 16 KDa

Observed band size: 16 KDa