
H2AFB1 Polyclonal Antibody

(Catalog # A62726)

Background

Atypical histone H2A which can replace conventional H2A in some nucleosomes and is associated with active transcription and mRNA processing. 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. Nucleosomes containing this histone are less rigid and organize less DNA than canonical nucleosomes in vivo. They are enriched in actively transcribed genes and associate with the elongating form of RNA polymerase. They associate with spliceosome components and are required for mRNA splicing. May participate in spermatogenesis.

Description

H2AFB1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P0C5Y9

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Histone H2A-Bbd type 1 protein (1-115AA)

Storage

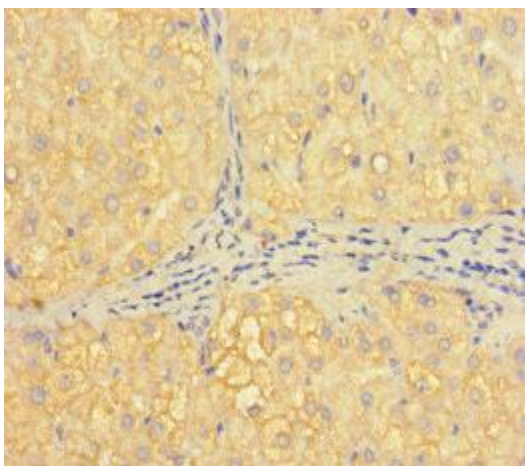
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

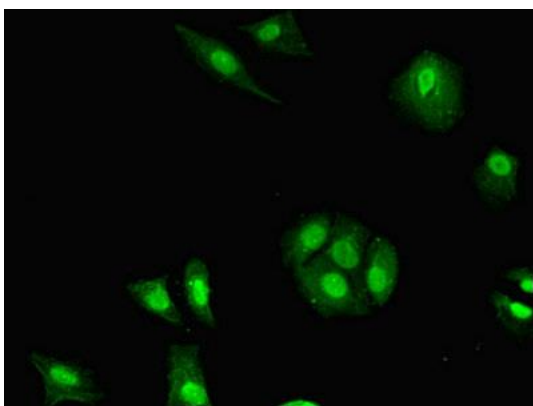
Histone H2A-Bbd type 1, H2A Barr body-deficient, H2A.Bbd, H2AFB1

Application

ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human liver tissue using H2AFB1 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of MCF-7 cells using H2AFB1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)