
Histone H2B Type 1-B Polyclonal Antibody

(Catalog # A50289)

Background

Core component of nucleosome. 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

Histone H2B Type 1-B Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P33778

Purification

Protein G purified

Immunogen

Recombinant Human Histone H2B type 1-B protein (2-126AA)

Storage

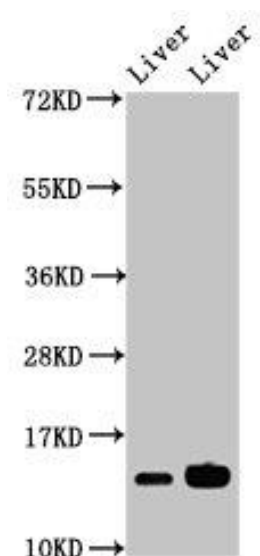
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Histone H2B.1, Histone H2B.f, H2B/f, H2BFF, HIST1H2BB

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:500-1:1000, IF:1:50-1:500



Western Blot

Positive WB detected in: Mouse liver tissue, Rat liver tissue

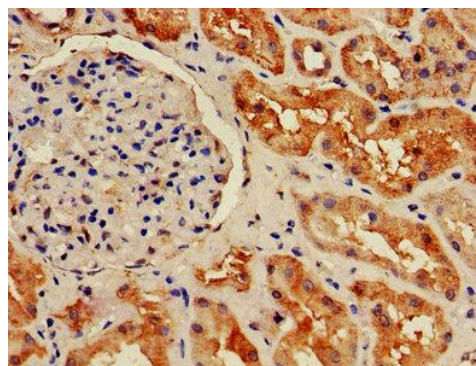
All lanes: HIST1H2BB antibody at 2.8ug/ml

Secondary

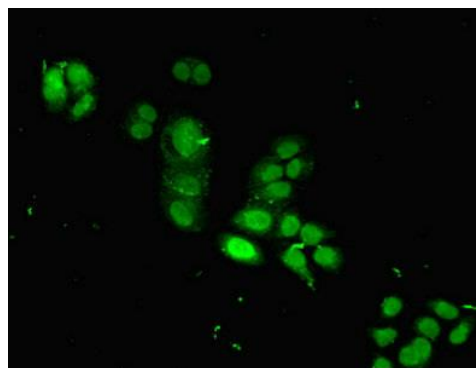
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 14 kDa

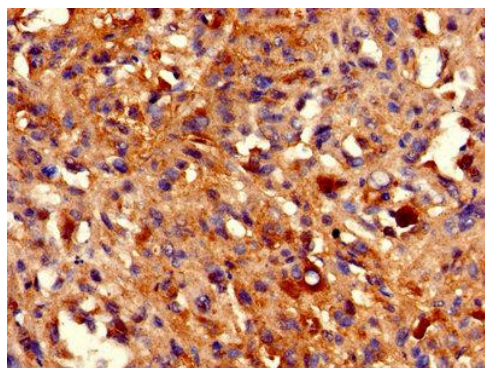
Observed band size: 14 kDa



IHC image of HIST1H2BB Antibody diluted at 1:800 and staining in paraffin-embedded human kidney tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Immunofluorescent analysis of PC-3 cells using HIST1H2BB Antibody at dilution of 1:266 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



IHC image of HIST1H2BB Antibody diluted at 1:800 and staining in paraffin-embedded human melanoma performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.