
Histone H2A.J Polyclonal Antibody

(Catalog # A53292)

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 H2A.J Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Preservative: 0.1mol/L NaCl, 15mmol/L NaN₃, PH7.2.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9BTM1

Purification

Protein G purified

Immunogen

Recombinant Human Histone H2A.J protein (2-129AA)

Storage

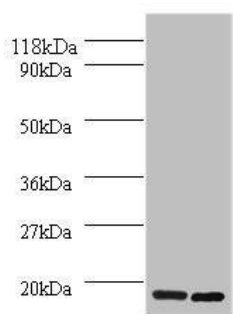
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

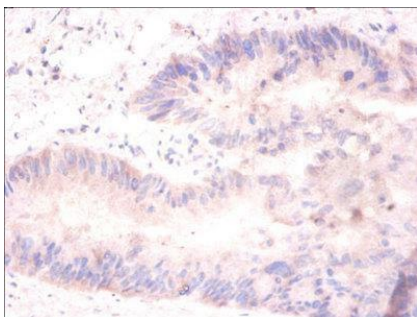
H2AFJ

Application

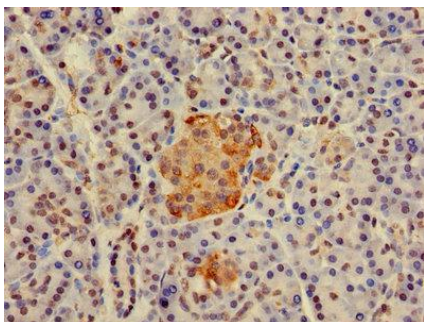
ELISA, WB, IHC; Recommended dilution: WB: 1:1000-1:5000, IHC:1:20-1:200



Western blot All lanes: H2AFJ antibody
at 2ug/ml Lane 1: Mouse brain tissue
Lane 2: MCF-7 whole cell lysate
Secondary
Goat polyclonal to rabbit IgG at 1/10000 dilution Predicted
band size: 15, 17 kDa
Observed band size: 15 kDa



Immunohistochemistry of paraffin-embedded human
colon cancer using H2AFJ Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human
pancreatic tissue using H2AFJ Antibody at dilution of
1:100