
Histone H2AX Polyclonal Antibody

(Catalog # A71359)

Background

Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. Two molecules of each of the four core histones (H2A, H2B, H3, and H4) form an octamer, around which approximately 146 bp of DNA is wrapped in repeating units, called nucleosomes. The linker histone, H1, interacts with linker DNA between nucleosomes and functions in the compaction of chromatin into higher order structures. This gene encodes a replication-independent histone that is a member of the histone H2A family, and generates two transcripts through the use of the conserved stem-loop termination motif, and the polyA addition motif.

Description

Histone H2AX Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat, Other (Wide Range)

Isotype

IgG

Uniprot ID

P16104

Purification

Affinity Purified

Immunogen

A synthetic phosphorylated peptide around S2 of human Histone H2AX (NP_002096.1)

Storage

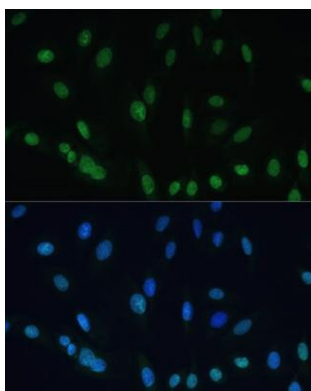
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

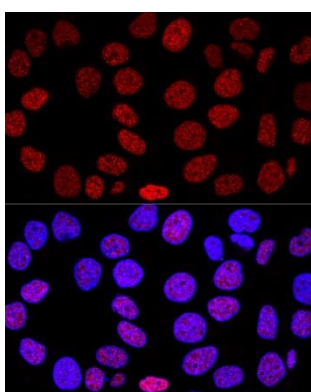
H2AFX; H2A.X; H2A/X; H2AX; histone H2AX

Application

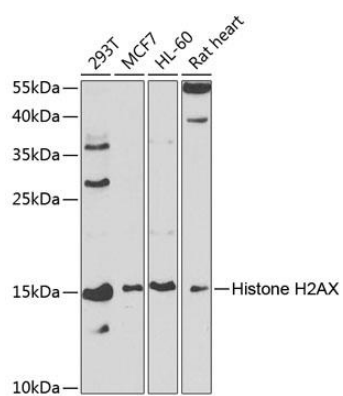
WB, IHC, IF; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:100, IF 1:50 - 1:100



Immunofluorescence analysis of U-2 OS cells using Histone H2AX Polyclonal Antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.



Confocal immunofluorescence analysis of HeLa cells using Histone H2AX Polyclonal Antibody at dilution of 1:200. Blue: DAPI for nuclear staining.



Western blot analysis of extracts of various cell lines, using Histone H2AX Polyclonal Antibody at 1:1000 dilution.

Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

Exposure time: 90s.