

Phospho Histone H2A (T121) Polyclonal Antibody

(Catalog #A68372)

Description

Phospho Histone H2A (T121) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

PBS with 50% glycerol, 0.5% BSA* and 0.02% sodium azide.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P0C0S8/P04908/P20671/Q96KK5/Q6FI13/Q7L7L0

Purification

Affinity Purified

Immunogen

Synthetic Peptide

Storage

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

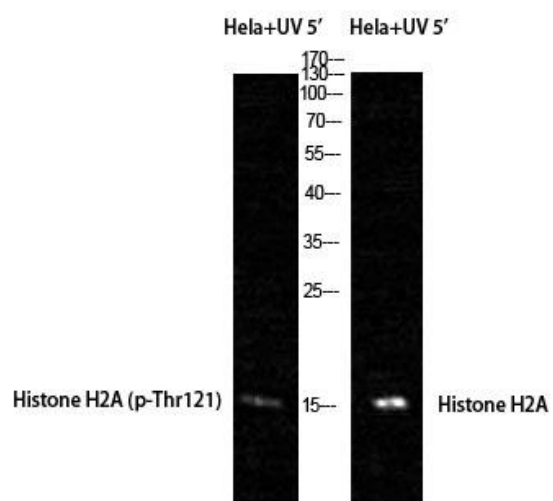
Alternative Names

HIST1H1A; H1F1; Histone H1.1; Histone H1a; HIST1H1C; H1F2; Histone H1.2; Histone H1c; Histone H1d; Histone H1s-1; HIST1H1D; H1F3; Histone H1.3; Histone H1c; Histone H1s-2; HIST1H1E; H1F4; Histone H1.4; Histone H1b; Histone H1s-4; H2AT121; H2AT121ph

Application

WB, IHC, ELISA; Recommended dilution: IHC:1:100-1:300, ELISA:1:5000

**The BSA is derived from animal or animal-derived material of negligible amounts. The animal-derived material is subject to heat treatment at a temperature higher than 65°C for at least three hours and acid treatment with pH value of less than 5 for at least three hours, thereby being free of Bovine Spongiform Encephalopathy (BSE) and Transmissible Spongiform Encephalopathy (TSE). Additionally, as all BSA-containing products are strictly intended for research purposes and not for diagnostic or therapeutic use, they are not subject to certification authority oversight.*



Western blot analysis of HeLa+UV 5' cells using the Phospho Histone H2A (T121) Polyclonal Antibody.