

Phospho Histone H2A.X (Tyr142) Polyclonal Antibody

(Catalog # A68363)

Description

Phospho Histone H2A.X (Tyr142) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

PBS with 0.02% sodium azide, 50% glycerol, pH 7.4.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P16104

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

H2AFX; H2AX; Histone H2A.x; H2a/x; H2A.XTyr142; H2A.XTyr142ph; H2A.XT142ph

Application

WB; Recommended dilution: WB: 1:1000-2000



Western blot analysis of extracts from HeLa cells, 1:2000.