

Histone H2A.ZK7ac (Acetyl H2A.ZK7) Polyclonal Antibody

(Catalog #A68351)

Description

Histone H2A.ZK7ac (Acetyl H2A.ZK7) 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

P0C0S5

Purification

Affinity Purified

Immunogen

Synthetic Peptide

Storage

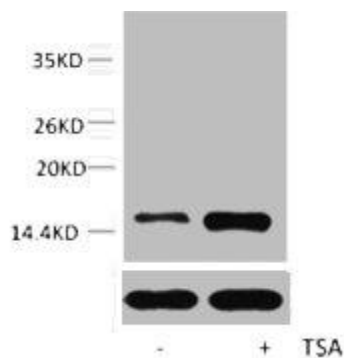
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

H2AFZ; H2AZ; Histone H2A.Z; H2A/z

Application

WB, ELISA; Recommended dilution: WB 1:500 - 1:1000, ELISA 1:20000



Western blot analysis of extracts from HeLa cells, untreated (-) or treated, 1:5000.