

Histone H2AS129ph antibody (pAb)

Catalog Nos: 39271, 39272

RRID: AB_2793213

Isotype: Serum

Application(s): ChIP, DB, WB

Reactivity: Budding Yeast, Wide Range Predicted

Volumes: 200 μ l, 10 μ l

Purification: None

Host: Rabbit

Molecular Weight: 14 kDa

Background: Histone H2A is one of the core components of the nucleosome, the basic building block of chromatin. Unlike mammals that have a distinct DNA damage-responsive histone H2A variant, H2AX, budding yeast histone H2A contains an SQ motif in the C-terminus that is phosphorylated by the Mec1 and Tel1 kinases in response to DNA double strand breaks. As with mammalian H2AX phosphorylation, phosphorylation of yeast histone H2A at serine 129 is required for the efficient repair of DNA double-strand breaks.

Immunogen: This Histone H2A phospho Ser129 antibody was raised against a peptide including phospho-serine 129 of budding yeast Histone H2A.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

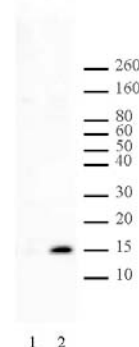
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:2,000 dilution

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

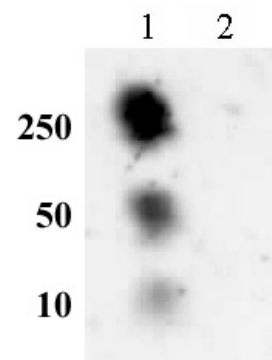


Histone H2AS129ph antibody (pAb) tested by Western blot.

Whole-cell extract of yeast exposed to 200Gy ionizing radiation blotted with Histone H2A phospho Ser129 pAb at a dilution of 1:1,000.

Lane 1: Yeast containing an alanine at position 129.

Lane 2: Wild-type yeast.



Histone H2AS129ph antibody (pAb) tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of Histone H2A phospho Ser129 pAb for phospho Ser129 histone H2A. Peptides corresponding to the immunogen and the unmodified version of the immunogen were spotted onto PVDF and probed with the antibody at 1:500. The amount of peptide (picomoles) spotted is indicated next to each row.

Lane 1: Phospho-Ser129 peptide.