

Htz1 / Histone H2A.Z antibody (pAb)

Catalog Nos: 39647, 39648

RRID: AB_2793289

Isotype: IgG

Application(s): ChIP, ChIP-Seq, WB

Reactivity: Budding Yeast

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 15 kDa

Background: Htz1 is the *S. cerevisiae* homolog of histone H2AZ and is involved in transcriptional regulation and chromosome architecture.

Immunogen: This Htz1 antibody was raised against full-length Htz1 protein from *S. cerevisiae*.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

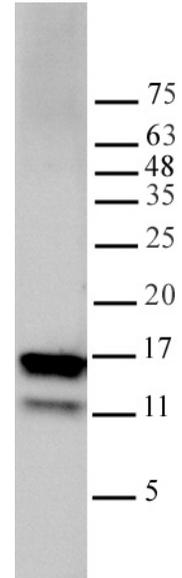
Application Notes:

Applications Validated by Active Motif:

WB: 1:300 - 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.



Htz1 pAb tested by Western blot.

Whole-cell extract from *S. cerevisiae* was probed with Htz1 / Histone H2A.Z antibody (pAb) at a 1:300