

Hbo1 antibody (pAb)

Catalog Nos: 39105, 39106

RRID: AB_2793156

Isotype: Serum

Application(s): WB

Reactivity: Human, Wide Range Predicted

Volumes: 200 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 75 kDa

Background: Hbo1 (HAT binding to ORC1) is a MYST-family histone acetyltransferase (HAT) that forms a complex with ING4/5, JADE1/2/3, and Eaf6, which is found enriched on coding regions of genes, supporting a role in transcriptional elongation. In addition, Hbo1 is thought to participate in events controlling the initiation of DNA replication.

Immunogen: This Hbo1 antibody was raised against a synthetic peptide containing amino acids 158-172 of human Hbo1.

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

Application Notes:

Applications Validated by Active Motif:

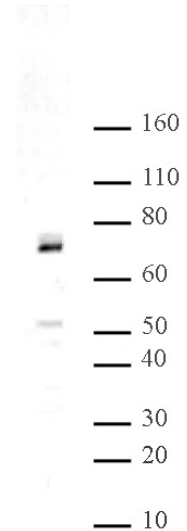
WB*: 1:500 - 1:2,000 dilution

For optimal results, primary antibody incubations should be performed overnight at 4°C. Individual optimization may be required.

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts.

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.



Hbo1 pAb tested by Western blot.

Detection of Hbo1 by Western blot analysis. HeLa whole-cell extract (15 µg) was probed with Hbo1 pAb at a 1:500 dilution.