

## LAP2 alpha antibody (mAb)

**Catalog Nos:** 39267, 39268

**RRID:** AB\_2793212

**Clone:** 3A3

**Isotype:** IgG1

**Application(s):** WB

**Reactivity:** Human

**Volumes:** 200 µl, 10 µl

**Purification:** Ascites

**Host:** Mouse

**Molecular Weight:** 80 kDa

**Background:** LAP2 α (Lamin Associated Polypeptide, alpha isoform, Thymopoietin) is a nucleoplasmic protein that oligomerizes and binds Hsp70 and Hsc70. It is involved in the organization of the nuclear structure through its interaction with A-type Lamins (Lamin B1). LAP2 α regulates the cell cycle through its interaction with the Rb protein.

**Immunogen:** This LAP2 alpha antibody was raised against a peptide including the C-terminus of human LAP2 alpha.

**Buffer:** Ascites containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

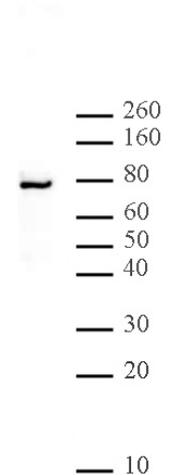
**Application Notes:**

Applications Validated by Active Motif:

WB: 1:1,000 - 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.



**LAP2 alpha mAb (Clone 3A3) tested by Western blot.**

Detection of LAP2 alpha by Western blot. The analysis was performed using HeLa nuclear extract and LAP2 alpha mAb (Clone 3A3) at a 1:2,000 dilution.