

## L3MBTL2 antibody (pAb)

**Catalog Nos:** 39569, 39570

**RRID:** AB\_2615062

**Isotype:** Serum

**Application(s):** IP, WB

**Reactivity:** Human

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** None

**Host:** Rabbit

**Molecular Weight:** 90 kDa

**Background:** L3MBTL2 (Lethal (3) malignant brain tumor-like 2 protein) is a putative member of the polycomb group family and is the human homolog of *Drosophila* Sfmtb, critical for Hox gene silencing. L3MBTL2 contains one FCS-type zinc finger and four MBT (malignant brain tumor) repeats. L3MBTL2 associates with chromatin-remodeling complexes and helps inhibit the expression of proteins that trigger the cells into mitosis. L3MBTL2 is a part of a complex involved in the G<sub>0</sub> phase of the cell cycle. This complex contains other proteins such as E2F-6, Max or HP1 gamma, and participates in transcriptional repression.

**Immunogen:** This L3MBTL2 antibody was raised against a His-Tagged fusion protein containing amino acids 1-230 of human L3MBTL2.

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

### Application Notes:

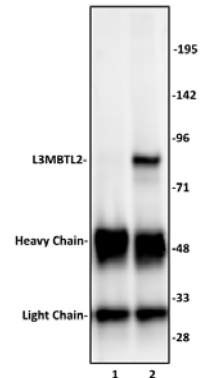
Applications Validated by Active Motif:

IP: 10  $\mu$ l per IP

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

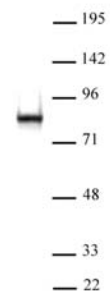
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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



### L3MBTL2 antibody (pAb) tested by Immunoprecipitation.

10  $\mu$ l of L3MBTL2 antibody was used to immunoprecipitate L3MBTL2 from 500  $\mu$ g of HeLa nuclear cell extract (lane 2). 10  $\mu$ l of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the L3MBTL2 antibody at a dilution of 1:2,000.



### L3MBTL2 pAb tested by Western blot.

Detection of L3MBTL2 by Western blot analysis with 20  $\mu$ g HeLa nuclear extract and a 1:2,000 dilution of L3MBTL2 pAb.