

| PRODUCT NAME                            |                         |                                 |
|---|-------------------------|---------------------------------|
| L3MBTL1 polyclonal antibody             |                         |                                 |
| Other names: L3MBTL, L3MBT              |                         |                                 |
| <b>Cat. No.</b> C15410023 (pAb-023-050) | <b>Type:</b> Polyclonal | <b>Size:</b> 50 µg/ 25 µl       |
| <b>Lot #:</b> 001                       | <b>Source:</b> Rabbit   | <b>Concentration:</b> 2.0 µg/µl |

**Description:** Polyclonal antibody raised in rabbit against human L3MBTL1 [lethal 3 malignant brain tumor-like protein], using the N-terminal part of the protein.

**Specificity:** Human: positive  
Other species: not tested

| Applications        | Suggested dilution | References |
|---------------------|--------------------|------------|
| Western blotting    | 1:1,000            | Fig1       |
| Immunoprecipitation | 1 - 5 µg per IP    | Fig 2      |

**Purity:** Protein G purified polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

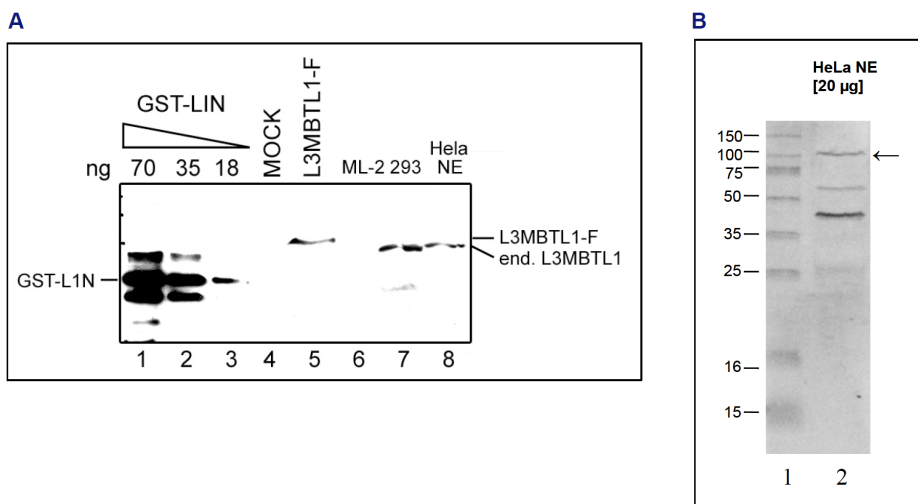
**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Last data sheet update:** April 9, 2010

#### Target description

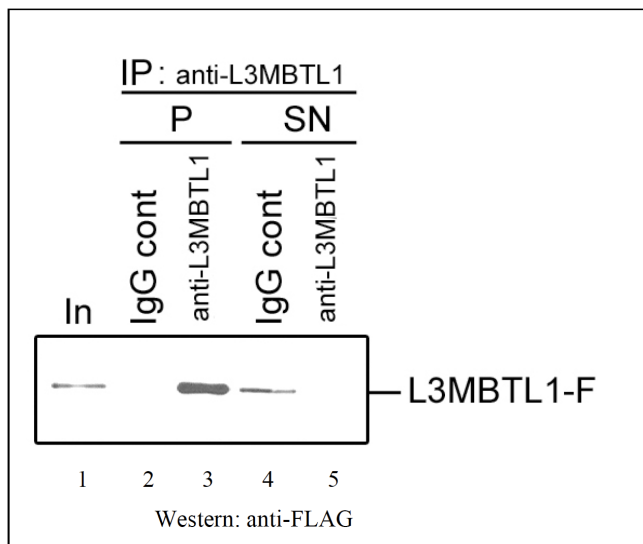
L3MBTL1 (UniProtKB/Swiss-Prot entry Q9Y468) is a member of the polycomb group (PcG) proteins, which function as transcriptional regulators in large protein complexes. In Drosophila, homozygous mutations in the L3MBTL1 gene have been demonstrated to cause brain tumors, identifying L3MBTL1 as a tumor suppressor gene. L3MBTL1 forms a homodimer and interacts with TEL (ETV6). L3MBTL1 probably plays a role in cell proliferation. Overexpression of L3MBTL1 induces multinucleated cells, suggesting that it is required for normal mitosis.



**Figure 1**

**Western blot analysis using the Diagenode antibody directed against L3MBTL1**

Western blot was performed on nuclear extracts from HeLa cells (HeLa NE) using the Diagenode antibody against L3MBTL1 (Cat. No. pAb-023-050) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk (lane 2). A molecular weight marker is shown in lane 1; the location of the protein of interest is indicated on the right.



**Figure 2**

**Immunoprecipitation using the Diagenode antibody directed against L3MBTL1**

Immunoprecipitation was performed on nuclear extracts from cells transfected with FLAG-tagged L3MBTL1 (L3MBTL1-F) with the Diagenode antibody against L3MBTL1 (Cat. No. pAb-023-050) and with an IgG negative control antibody. Western Blot was performed with an anti-FLAG antibody to demonstrate the presence of L3MBTL1-F in the input (lane 1), in the precipitated fraction (P) and in the supernatant (SN). These data show that L3MBTL1-F was efficiently precipitated by the L3MBTL1 antibody (lane 3 and 5), whereas it was not precipitated by the IgG negative control antibody (lane 2 and 4).