

# Anti-ILPIP (IN)

## CATALOG NO.: 54150

## BACKGROUND:

ILPIP (ILP-interacting protein) has recently been shown to interact with XIAP, a member of the IAP (Inhibitor of Apoptosis) protein family (1). Together, these two proteins synergistically activate the TAK1/JNK1 signal transduction pathway, which acts to protect against the interleukin-1β converting enzyme of Fas-induced apotosis (1,2). This protein was independently isolated and characterized as polyploidy-associated protein kinase (PAPK), a member of the Ste20/germinal center kinase family that modulates cytoskeletal organization and cell survival (3), suggesting that its protective role may be a function of its kinase activity. Although two mRNA sequences have been isolated, only the shorter isoform has been observed.

#### SOURCE & REACTIVITY:

Rabbit anti-ILPIP polyclonal antibody was raised against a synthetic peptide corresponding to amino acids near the center of human ILPIP (Genbank accession NP\_061041). Anti-ILPIP reacts with ILPIP at the molecular weight of 34 kDa on western blot. Species reactivity includes human, mouse, and rat, while others are not tested.

## **APPLICATION:**

The following concentration ranges are recommended starting points for this product.



Positive Control: Human brain cell lysate



Immunohistochemistry of ILPIP in human brain tissue with anti-ILPIP at 10 µg/ml.

This product is for in vitro research purposes only.

#### **RELATED PRODUCTS:**

Anti-ILPIP (CT), Catalog No. **54151** Anti-XIAP (CT), Catalog No. **54174** Anti-cIAP (CT), Catalog No. **54173** 

## STORAGE:

The antibody is supplied as immunoaffinity purified IgG in 1X PBS containing 0.02% sodium azide. Store at 4°C, stable for one year. Store at 2-8 °C for up to 1 year. Avoid repeated freeze thaw cycles.

#### **REFERENCES:**

- 1. Sanna MG. et al. J. Biol. Chem. 277, 30454(2002).
- 2. Sanna MG. et al. Proc. Natl. Acad. Sci. USA 95 6015(1998).
- 3. Nishigaki K. et al. J. Biol. Chem. 278, 13520(2003).