

Anti-Toso

(Regulator of Fas-induced apoptosis, Fas apoptotic inhibitory molecule 3, FAIM3)

CATALOG NO.: 54102

BACKGROUND:

Apoptosis is an important process by which normal tissue homeostasis and function are maintained. One of the major signals that regulate this process is mediated by the activation of the Fas receptor by its ligand. This leads to the formation of a Fas-associated death domain (FADD)- containing death-inducing signaling complex and the activation of caspase-8, which in turn activates downstream effector caspases, such as caspase-3 and -7 (1). Recent experiments have shown that over-expression of Toso, a novel regulator of Fas-induced apoptosis in lymphoid cells, in Jurkat cells as well as transgenic mice render these cells resistant to Fas-induced apoptosis, but not to TRAIL-induced apoptosis (2,3). Furthermore, Toso was found to associate with FADD, suggesting that Toso functions by disrupting the formation of the death-inducing signaling complex.

SOURCE & REACTIVITY:

Rabbit anti-Toso polyclonal antibody was raised against a 13 amino acid peptide from near the C-terminus of human Toso (GenBank accession no. NP_005440). Anti-Toso reacts with Toso at the molecular weight of 60 kDa on western blot. Species reactivity includes human, while others are not tested.

APPLICATION:

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

WB: 1 µg/ml.

Positive Control: Human lung tissue lysate



This product is for in vitro research purposes only.

RELATED PRODUCTS:

Anti-TRAIL (CT), Catalog No. **54008** Anti-Caspase-7 (CT), Catalog No. **54211** Anti-Caspase-8 (CT), Catalog No. **54214**

STORAGE:

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

REFERENCES:

- 1. Curtin, JF. et al. Cell Signal. 15, 983 (2003).
- 2. Hitoshi, Y. et al. Immunity 8, 461 (1998).
- 3. Song, Y. et al. J. Biol. Chem. 280, 9618 (2005).