

Anti-XAF-1 (CT) (XIAP associated)

CATALOG NO.: 54153

BACKGROUND:

XAF-1 binds to XIAP, an inhibitor of caspases-3, -7, and – 9, and triggers its relocation from the cytosol to the nucleus (1, 2). Over-expression of XAF-1 results in the neutralization of XIAP's ability to inhibit cell death (1). XAF-1 is normally expressed in all adult and fetal tissues, but was found to be present in very low levels in a variety of cancer cell lines (3). In contrast, XIAP levels have been shown to be high in a majority of cell lines. Low XAF-1 and high basal expression of XIAP may play a critical role in maintaining survival of cancer cell lines (2, 3). Both IFN- α 2 and IFN- β can induce XAF-1 mRNA in all cells examined but induction of XAF-1 protein (as observed by immunoblot analysis) was seen only in cell lines sensitive to the apoptotic effects of IFNs (4).

SOURCE & REACTIVITY:

Rabbit anti-XAF-1 polyclonal antibody was raised against a synthetic peptide corresponding to amino acids at the Cterminus of human XAF-1 (Genbank accession CAA68030). Anti-XAF-1 reacts with XAF-1 at the molecular weight of 32 kDa on western blot. Species reactivity includes human and mouse, while others are not tested.

APPLICATION:

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

WB: 0.5 to 2 µg/ml.

Positive Control: Human spleen cell lysate



Immunohistochemical staining of human spleen tissue using anti-XAF-1 at 2 μ g/ml.



This product is for in vitro research purposes only.

RELATED PRODUCTS:

Anti-XAF-1 (IN), Catalog No. **54192** Anti-XIAP (CT), Catalog No. **54174**

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. Liston, P. et al. Nature Cell Biol. 3, 128 (2001).
- 2. Deveraux, QL. et al. Nature 388, 300 (1997).
- 3. Fong, WG. et al. Genomics 70, 113 (2000).
- 4. Leaman, DW. et al. J. Biol. Chem. 277, 28504 (2002).