

## Anti-XIAP

Rabbit Polyclonal Antibody

**Catalog # X08-63R**

Lot # O2109-24

### Cited Applications

WB, ELISA, IF, IHC

*Ideal working dilutions for each application should be empirically determined by the investigator.*

### Specificity

Recognizes the XIAP protein

### Cross Reactivity

Human and Mouse

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

XIAP antibody was raised against a synthetic peptide corresponding to 13 amino acids at the C-terminus of human XIAP

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

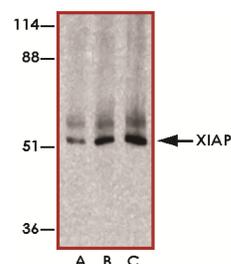
### Scientific Background

XIAP is a member of a family of proteins which inhibit apoptosis and act as mammalian cell-death suppressors. XIAP inhibits at least 2 members of the caspase family of cell-death proteases, caspase-3 and caspase-7 (1). In addition, XIAP inhibits apoptosis induced by menadione, a potent inducer of free radicals, and ICE. Furthermore, XIAP can mediate protection of cells from apoptosis by utilizing both a JNK1 activation pathway that involves ILPIP and a caspase inhibition pathway that is independent of ILPIP (2). Disruption of the XIAP gene in human colon cancer cells can enhance sensitivity of these cells to exogenously added TRAIL suggesting that XIAP is a nonredundant modulator of TRAIL-mediated apoptosis.

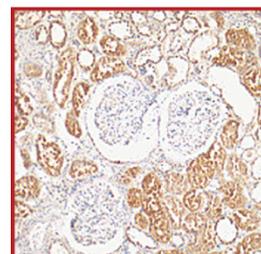
### References

1. Deveraux Q L, et al: X-linked X-linked IAP is a direct inhibitor of cell-death proteases. Nature 388: 300-304, 1997.
2. Tang G, et al: Inhibition of JNK activation through NF-kappa-B target genes. Nature 414: 313-317, 2001.

### Sample Data



Western blot analysis of XIAP in human kidney lysate with XIAP antibody at 0.5 (lane A), 1 (lane B), and 2 (lane C) µg/ml, respectively.



Immunohistochemical staining of human kidney tissue using XIAP antibody at 2 µg/ml.

## Anti-XIAP

Rabbit Polyclonal Antibody

Catalog Number

X08-63R

Specific Lot Number

O2109-24

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**