

## Anti-BAX

Rabbit Polyclonal Antibody

**Catalog # B14-363R**

Lot # O2109-28

### Cited Applications

WB, ELISA, ICC, IF

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

### Specificity

Recognizes the BAX protein

### Cross Reactivity

Human

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

BAX antibody was raised against a peptide corresponding to 16 amino acids near the amino-terminus of human BAX

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

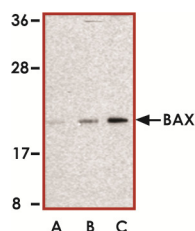
### Scientific Background

BAX is a proapoptotic protein of the BCL2 protein family. BAX forms a heterodimer with BCL2 and functions as an apoptotic activator. BAX interacts with and increases the opening of the mitochondrial voltage-dependent anion channel (VDAC), which leads to the loss in the mitochondrial membrane potential and the release of cytochrome c (1). The expression of BAX gene is regulated by the tumor suppressor p53 and BAX has been shown to be involved in p53-mediated apoptosis. Multiple alternatively spliced transcript variants, which encode different isoforms, have been reported for this gene (2).

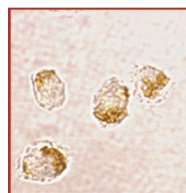
### References

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 BAX in HL-60 cell lysates with BAX antibody at (A) 1, (B) 2, and (C) 4 µg/ml.



Immunocytochemistry staining of HL-60 cells using BAX at 2 µg/ml.

## Anti-BAX

Rabbit Polyclonal Antibody

Catalog Number

**B14-363R**

Specific Lot Number

**O2109-28**

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.**