

Anti-BCL2

Rabbit Polyclonal Antibody

Catalog # B25-63CR

Lot # O2109-26

Cited Applications

WB, ELISA, ICC

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

Specificity

Recognizes the BCL2 protein

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

BCL2 antibody was raised against a peptide corresponding to 14 amino acids near the middle of human BCL2

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

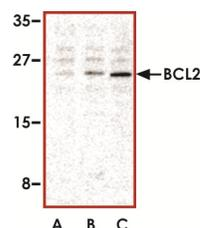
Scientific Background

BCL2 gene encodes an integral inner mitochondrial membrane protein that acts as an antiapoptotic protein (1). The protein BAD can antagonize both the cell cycle and antiapoptotic functions of BCL2 through binding to the BH3 domain. Constitutive expression of BCL2, such as in the case of translocation of BCL2 to Ig heavy chain locus, is thought to be the cause of follicular lymphoma (2). BCL2 is phosphorylated on specific serine/threonine residues within the unstructured loop in response to diverse stimuli and such phosphorylation has been associated with the loss of the biological function of BCL2.

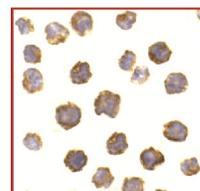
References

1. Hockenbery, D. et al: Bcl-2 is an inner mitochondrial membrane protein that blocks programmed cell death. *Nature* 348: 334-336, 1990.
2. Tsujimoto, Y. et al: Involvement of the bcl-2 gene in human follicular lymphoma. *Science* 228: 1440-1443, 1985.

Sample Data



Western blot analysis of BCL2 in A-20 cell lysates with BCL2 antibody at (A) 1, (B) 2, and (C) 4 µg/ml.



Immunocytochemistry of BCL2 in A20 cells with BCL2 antibody at 2 µg/ml.

Anti-BCL2

Rabbit Polyclonal Antibody

Catalog Number

B25-63CR

Specific Lot Number

O2109-26

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
www.signalchem.com

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