

Anti-BCL2

Rabbit Polyclonal Antibody

Catalog #**B25-63R**

Lot # H245-7

Cited Applications

- Immunoprecipitation (2 µg)
- Western blot (1:1000)

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

Specificity

Recognizes the BCL2 protein

Cross Reactivity

- Immunoprecipitation of BCL2 from human cells
- Western blot of BCL2 from human cells

BCL2 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 54-58 of human BCL2

Formulation

PBS pH 7.4, 0.02% sodium azide, 50% glycerol

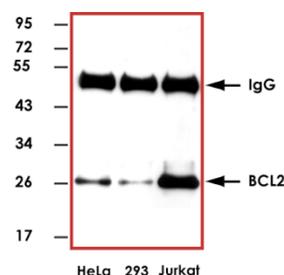
Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. 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.

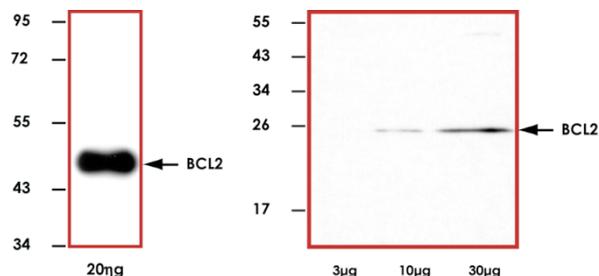
Scientific Background

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.

Sample Data



Representative immunoprecipitation with 2 µg of Anti-BCL2 using 500 µg of HeLa, HEK-293T, and Jurkat cell lysate.



Representative western blot with Anti-BCL2 (1:1000) using 20 ng of GST tagged human recombinant BCL2 (left image) or 3 µg, 10 µg, and 30 µg of Jurkat cell lysate.

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.

Anti-BCL2

Rabbit Polyclonal Antibody

Catalog Number	B25-63R
Specific Lot Number	H245-7
Purification	Affinity chromatography
Concentration	1.0 µg/µL
Stability	1 yr 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.