

Anti-BCL2L14

Rabbit Polyclonal Antibody

Catalog # B26-363NR

Lot # O2361-16

Cited Applications

E, WB, IHC, IF

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

Specificity

Recognizes the BCL2L14 protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

BCL2L14 antibody was raised against a peptide corresponding to 15 amino acids near the C-terminus of human BCL2L14

Formulation

PBS + 0.02% sodium azide.

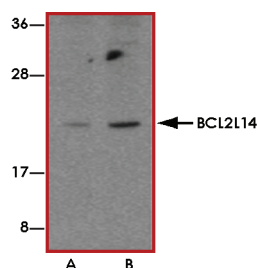
Stability

1yr at -20°C from date of shipment.

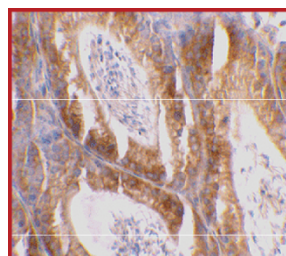
References

1. Cory S et al: The Bcl-2 family: roles in cell survival and oncogenesis. *Oncogene* 2003; 22:8590-607
2. Heiser D et al: The Bcl-2 protein family and its role in the development of neoplastic disease. *Exp. Geront.* 2004; 39:1125-35.
3. Guo B et al: Bcl-G, a novel pro-apoptotic member of the Bcl-2 family. *J. Biol. Chem.* 2000; 276:2780-5.

Sample Data



Western blot analysis of BCL2L14 in U937 cell lysates with BCL2L14 antibody at (A) 2.5 and (B) 5 µg/mL.



Immunohistochemical staining of mouse testis tissue using BCL2L14 antibody at 2 µg/mL.

Anti-BCL2L14

Rabbit Polyclonal Antibody

Catalog Number

B26-363NR

Specific Lot Number

O2361-16

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.