
Bcl2 Polyclonal Antibody

(Catalog #A52548)

Background

Suppresses apoptosis in a variety of cell systems including factor-dependent lymphohematopoietic and neural cells. Regulates cell death by controlling the mitochondrial membrane permeability. Appears to function in a feedback loop system with caspases. Inhibits caspase activity either by preventing the release of cytochrome c from the mitochondria and/or by binding to the apoptosis-activating factor (APAF-1).

Description

Bcl2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300. 50% Glycerol, 0.01M PBS, pH 7.4.

Specificity

Mouse

Isotype

IgG

Uniprot ID

P10417

Purification

Protein G purified

Immunogen

Recombinant Mouse Apoptosis regulator Bcl-2 protein (5-205AA)

Storage

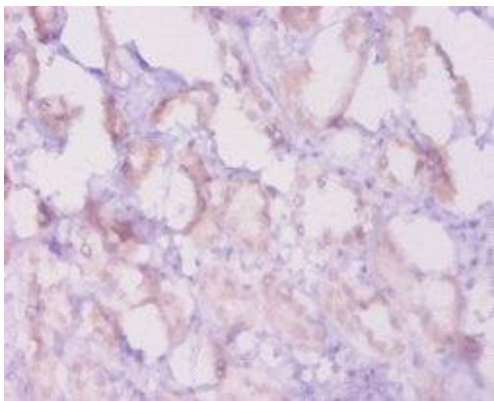
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Bcl2, Bcl-2

Application

ELISA, IHC; Recommended dilution: IHC:1:10-1:100



Immunohistochemistry of paraffin-embedded human kidney at dilution of 1:20