

PE Mouse Anti-human BCL2 Antibody
Bcl-2/100, monoclonal, Cross AdsorbedCatalog number: V103160
Unit size: 100 tests**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	BCL2
Clone	Bcl-2/100
Conjugate	PE

Biological Properties

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse serum and then conjugated with PE under optimal conditions
Application	FC (QC TESTED)

Applications

Apoptosis regulator Bcl-2 is a 26 kDa transmembrane protein that can be located in the cytosol, nuclear membrane and mitochondrion of cells. In Homo sapiens, apoptosis regulator Bcl-2 has been closely linked to critical functions such as protein heterodimerization activity. Apoptosis regulator Bcl-2 negatively regulates osteoblast proliferation, mitochondrial depolarization and cellular pH reduction. But it also positively regulates catalytic activity, B cell proliferation and cell population proliferation. It binds to repressing transcription factor, sequence-specific DNA and ubiquitin protein ligase. Apoptosis regulator Bcl-2 is the subject of extensive application due to the fact that it is a member of the negative regulation of intrinsic apoptotic signaling pathway, extrinsic apoptotic signaling pathway via death domain receptors and cytokine-mediated signaling pathway. Apoptosis regulator Bcl-2 takes part in processes such as organismal processes, for example, T cell differentiation in thymus, response to UV-B and ovarian follicle development. It is thought to be essential to glycoprotein biosynthetic process, mitochondrial membrane potential and calcium ion transport.