

BCL-2 (Ab-69) Antibody [#B0774]

Catalog Number:	B0774
Concentration:	1mg/ml
Swiss-Prot No.:	P10415
Other Names:	Apoptosis regulator Bcl-2, BCL2
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human BCL-2 around the phosphorylation site of threonine 69 (A-R-T ^P -S-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BCL-2 (Ab-69) antibody detects endogenous levels of total BCL-2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	H. Kurosawa, Am J Physiol Gastrointest Liver Physiol, Jun 1997; 272: 1587. JR Park, Blood, Aug 1995; 86: 868 - 876. A Plettenberg, Am. J. Pathol., Mar 1995; 146: 651. JM Ben-Ezra, Am. J. Pathol., Nov 1994; 145: 1036.