

BCL-2 (Ab-70) Antibody [#B7025]

Catalog Number:	B7025
Concentration:	1mg/ml
Swiss-Prot No.:	P10415
Other Names:	Apoptosis regulator Bcl-2; B-cell CLL/lymphoma 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 serine 70 (R-T-S ^P -P-L).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BCL-2 (Ab-70) antibody detects endogenous levels of total BCL-2 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	Tsujimoto Y., Proc. Natl. Acad. Sci. U.S.A. 83:5214-5218(1986). Eguchi Y., Nucleic Acids Res. 20:4187-4192(1992). Cleary M.L., Cell 47:19-28(1986).