

BCLW Antibody [#C13031]

Catalog Number:	C13031
Concentration:	1mg/ml
Swiss-Prot No.:	Q92843
Other Names:	Apoptosis regulator Bcl-W, Bcl-2-like 2 protein, BCL2L2, BCLW, KIAA0271
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 peptide derived from internal of human BCLW.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BCLW antibody detects endogenous levels of total BCLW protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
References:	Gibson L., Oncogene 13:665-675(1996). Nagase T., DNA Res. 3:321-329(1996). O'Reilly L.A., Cell Death Differ. 8:486-494(2001).