

BCL-XL (Ab-62) Antibody [#B7027]

- Catalog Number:** B7027-1, B7027-2, B7027
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** This antiserum was produced against synthesized non-phosphopeptide derived from human BCL-XL around the phosphorylation site of serine 62.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** BCL-XL (Ab-62) antibody detects endogenous levels of total BCL-XL protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:10000
- Swiss-Prot No. :** Q07817
- References:** Basu A, et.al.(2003) Mar 13; 538(1-3): 41-7.
Cheng E.H.-Y. et.al. (1996) Nature 379:554-556.
Sattler M. et.al. (1997), Science 275:983-986.
Muchmore S.W.et.al (1996).Nature 381:335-341.
Yu J., Zhang L. Mol. (2001)Cell 7:673-682.