



BCL-XL (Ab-62) Antibody

Catalog Number 21061

(For Research Use Only)

- Catalog Number:** 21061-1, 21061-2, 21061
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.

