

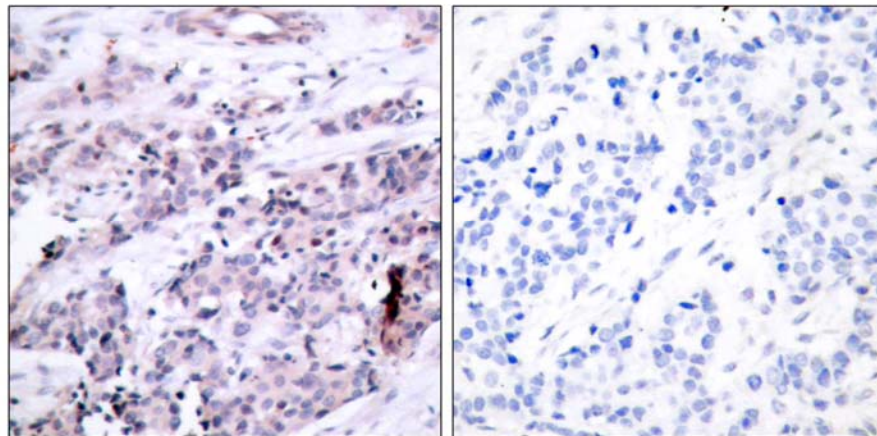


BCL-2(Ab-56) Antibody

Catalog Number 21059

(For Research Use Only)

- Catalog Number:** 21059-1, 21059-2, 21059
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-2 around the phosphorylation site of Threonine 56.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: BCL-2(Ab-56) Antibody detects endogenous levels of BCL-2 protein around Threonine 56.
Reactivity: Human
Applications: IHC: 1:50~1:100 ELISA: 1:10000
Swiss-Prot No. : P10415
References: Ling, Y. H. et al. (1998) *J. Biol. Chem.* 273, 18984-18991.
Huang, S.J. and Cidlowski, J.A. (2002) *FASEB* 16, 825-832.
Deng, X. et al. (2001) *J. Biol. Chem.* 276, 23681-23688.
Huang ST, et al. (2002) *FASEB J Jun*; 16(8): 825-32
Yamamoto, K. et al. (1999) *Mol. Cell. Biol.* 19, 8469-8478.



Immunohistochemical analysis of paraffin-embedded breast carcinoma, using BCL-2(Ab-56) Antibody (#21059). Left: Untreated; Right: Treated with synthesized peptide.