

14-3-3 ζ/δ (Ab-232) Antibody [#B0759]

Catalog Number:	B0759
Concentration:	1mg/ml
Swiss-Prot No.:	P63104
Other Names:	14-3-3 protein zeta/delta, 1433Z, 143Z, FAS, Factor activating exoenzyme S, KCIP-1, Mitochondrial import stimulation factor S1 subunit, Protein kinase C inhibitor protein-1, YWHAZ
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 14-3-3 ζ/δ around the phosphorylation site of threonine 232 (S-D-T ^P -Q-G).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	14-3-3 ζ/δ (Ab-232) antibody detects endogenous levels of total 14-3-3 ζ/δ protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Alastair Aitken, J. Biol. Chem., Mar 1995; 270: 5706. L Kockel, Genes & Dev., May 1997; 11: 1140 - 1147. Martijn J. van Hemert, J. Cell Sci., Mar 2004; 117: 1411 - 1420. Anne-Sophie Vercoutter-Edouart, Cancer Res., Jan 2001; 61: 76 - 80.