

## 14-3-3 $\zeta$ (Ab-58) Antibody [#B0001]

- Catalog Number:** B0001-1, B0001-2, B0001
- Amount:** 50 $\mu$ g/50 $\mu$ l, 100 $\mu$ g/100 $\mu$ l, 200 $\mu$ g/200 $\mu$ l
- Swiss-Prot No. :** P63104
- All 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
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg<sup>2+</sup> and Ca<sup>2+</sup>), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human 14-3-3  $\zeta$  around the phosphorylation site of serine 58 (R-S-S<sup>P</sup>-W-R).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** 14-3-3  $\zeta$  (Ab-58) antibody detects endogenous levels of total 14-3-3  $\zeta$  protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1000 IHC: 1:50~100 IP: Various Dilution  
IF: 1:500~1:1000 ELISA: 1:40000
- References:** Ying H. Shen Mol. Biol. Cell, Nov 2003; 14: 4721.  
David W. Powell, Mol. Cell. Biol., Aug 2003; 23: 5376 - 5387.  
Mackintosh. C, (2004) Biochem. J. 381, 329 - 42.  
Dougherty, M.K. and Morrison, D.K. (2004) J. Cell Sci. 117, 1875 - 84.