



## 14-3-3 $\zeta$ (Phospho-Ser58) Antibody

Catalog Number 11181

(For Research Use Only)

- Catalog Number:** 11181-1, 11181-2, 11181  
**Amount:** 50 $\mu$ g/50 $\mu$ l, 100 $\mu$ g/100 $\mu$ l, 200 $\mu$ g/200 $\mu$ l  
**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 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 phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.  
**Specificity/Sensitivity:** 14-3-3  $\zeta$  (Phospho-Ser58) antibody detects endogenous levels of 14-3-3  $\zeta$  only when phosphorylated at serine 58.  
**Reactivity:** Human, Mouse, Rat  
**Applications:** WB: 1:500~1:1000 IHC: 1:50 ~ 1:100  
**Swiss-Prot No. :** P10636  
**References:** Gu YM, et al. (2006) FEBS Lett ; 580(1): 305-310  
Powell DW, et al. (2003) Mol Cell Biol; 23(15): 5376-5387  
Mackintosh C. (2004) Biochem J; 381(Pt 2): 329-342.

