



14-3-3 ζ (Ab-58) Antibody

Catalog Number 21188

(For Research Use Only)

Catalog Number: 21188-1, 21188-2, 21188

Amount: 50 μ g/50 μ l, 100 μ g/100 μ l, 200 μ g/200 μ l

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 $^{\circ}$ C /1 year

Immunogen: The antiserum was produced against synthesized non-phosphopeptide derived from human 14-3-3 ζ around the phosphorylation site of serine 58 (R-S-S^P-W-R).

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.

Specificity/Sensitivity: 14-3-3 ζ (Ab-58) antibody detects endogenous levels of total 14-3-3 ζ protein.

Reactivity: Human, Mouse, Rat

Applications: WB: 1:500~1:1000 IHC: 1:50 ~ 1:100

Swiss-Prot No. : P63104

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.

