

## 14-3-3 $\beta$ Antibody [#C12001]

<b>Catalog Number:</b>	C12001
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P31946
<b>Other Names:</b>	14-3-3 protein beta/alpha, 1433B, 143B, KCIP-1, Protein 1054, Protein kinase C inhibitor protein-1, YWHAB
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	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.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human 14-3-3 $\beta$ .
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	14-3-3 $\beta$ antibody detects endogenous levels of total 14-3-3 $\beta$ protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Leffers H., J. Mol. Biol. 231:982-998(1993). Deloukas P., Nature 414:865-871(2001). Gevaert K., Nat. Biotechnol. 21:566-569(2003). Chi A., J. Proteome Res. 5:3135-3144(2006).