

14-3-3 θ Antibody [#C12006]

Catalog Number:	C12006
Concentration:	1mg/ml
Swiss-Prot No.:	P27348
Other Names:	14-3-3 protein T-cell, 14-3-3 protein tau, 1433T, 143T, HS1 protein, YWHAQ
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 peptide derived from internal of human 14-3-3 θ .
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	14-3-3 θ antibody detects endogenous levels of total 14-3-3 θ protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Nielsen P.J., Biochim. Biophys. Acta 1088:425-428(1991). Leffers H., J. Mol. Biol. 231:982-998(1993). Gevaert K., Nat. Biotechnol. 21:566-569(2003).