

14-3-3 θ/τ (Ab-232) Antibody [#B0758]

Catalog Number:	B0758
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 non-phosphopeptide derived from human 14-3-3 θ/τ around the phosphorylation site of serine 232 (S-D-S ^P -A-G).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	14-3-3 θ/τ (Ab-232) antibody detects endogenous levels of total 14-3-3 θ/τ protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Nielsen P.J., <i>im. Biophys. Acta</i> 1088:425-428(1991). Leffers H., <i>J. Mol. Biol.</i> 231:982-998(1993). Gevaert K., <i>Nat. Biotechnol.</i> 21:566-569(2003).