

14-3-3 η Antibody [#C12003]

Catalog Number:	C12003
Concentration:	1mg/ml
Swiss-Prot No.:	Q04917
Other Names:	14-3-3 protein eta, Protein AS1, YWHAH, YWHA1
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 η .
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	14-3-3 η antibody detects endogenous levels of total 14-3-3 η protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Swanson K.D., Biochim. Biophys. Acta 1216:145-148(1993). Ichimura-Ohshima Y., J. Neurosci. Res. 31:600-605(1992). Muratake T., Mol. Neurobiol. 11:223-230(1995).