

RIN1 Antibody [#C18175]

Catalog Number:	C18175
Concentration:	1mg/ml
Swiss-Prot No.:	Q13671
Other Names:	Ras and Rab interactor 1; Ras inhibitor JC99; Ras interaction/interference protein 1
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 RIN1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RIN1 antibody detects endogenous levels of total RIN1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Colicelli J., Proc. Natl. Acad. Sci. U.S.A. 88:2913-2917(1991). Han L., Mol. Cell. Biol. 15:1318-1323(1995). Han L., Proc. Natl. Acad. Sci. U.S.A. 94:4954-4959(1997).