

RHG22 Antibody [#C18371]

Catalog Number:	C18371
Concentration:	1mg/ml
Swiss-Prot No.:	Q7Z5H3
Other Names:	RHG22; Rho GTPase activating 2; Rho GTPase activating 22; RHOGAP2
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 RHG22.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RHG22 antibody detects endogenous levels of total RHG22 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Richnau N., J. Biol. Chem. 276:35060-35070(2001). Ota T., Nat. Genet. 36:40-45(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004).