

RHG17 Antibody [#C18367]

Catalog Number:	C18367
Concentration:	1mg/ml
Swiss-Prot No.:	Q68EM7
Other Names:	NADRIN; neuron-associated developmentally regulated protein; RHG17; rho GTPase activating protein 17; RhoGAP interacting with CIP4 homologs 1; RICH1; RICH1B; WBP15
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 RHG17.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RHG17 antibody detects endogenous levels of total RHG17 protein.
Reactivity:	Human, Mouse, Rat
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 German cDNA consortium; Submitted (JUL-2002) to the EMBL/GenBank/DDBJ databases.