

RGAP1 Antibody [#C11673]

Catalog Number:	C11673
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H0H5
Other Names:	Rac GTPase-activating protein 1, MgcRacGAP, RACGAP1, KIAA1478, MGCRCACGAP
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 RGAP1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RGAP1 antibody detects endogenous levels of total RGAP1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Kawashima T., Blood 96:2116-2124(2000). Wiemann S., Genome Res. 11:422-435(2001). Ota T., Nat. Genet. 36:40-45(2004).