

RAB3GAP1 Antibody [#C18124]

Catalog Number:	C18124
Concentration:	1mg/ml
Swiss-Prot No.:	Q15042
Other Names:	P130; RAB3 GTPase activating protein 1 (catalytic); RAB3 GTPase-activating protein 130 kDa; Rab3 GTPase-activating protein catalytic; Rab3-GAP; Rab3-GAP p130; RAB3GAP; RAB3GAP130; RB3GP; WARBM1
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 RAB3GAP1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RAB3GAP1 antibody detects endogenous levels of total RAB3GAP1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Nomura N., DNA Res. 1:223-229(1994). The German cDNA consortium; Submitted (SEP-2004) to the EMBL/GenBank/DDBJ databases. The MGC Project Team; Genome Res. 14:2121-2127(2004).