

RAB5C Antibody [#C18265]

Catalog Number:	C18265
Concentration:	1mg/ml
Swiss-Prot No.:	P51148
Other Names:	L1880; RAB5L; Ras-related protein Rab-5C
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 RAB5C.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RAB5C antibody detects endogenous levels of total RAB5C protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Han H.J., Cytogenet. Cell Genet. 73:137-139(1996). Albertsen H.M., Nat. Genet. 7:472-479(1994). Clemens D.L., Infect. Immun. 68:2671-2684(2000).