

RAP2C Antibody [#C18282]

Catalog Number:	C18282
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y3L5
Other Names:	RAP2C, member of RAS oncogene family; Ras-related protein Rap-2c
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 RAP2C.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RAP2C antibody detects endogenous levels of total RAP2C protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Shan Y.X., Submitted (MAY-2003) to the EMBL/GenBank/DDBJ databases. Ross M.T., Nature 434:325-337(2005). Paganini S., Biochimie 88:285-295(2006).