

RapGEF1 Antibody [#B1120]

Catalog Number:	B1120
Concentration:	1mg/ml
Swiss-Prot No.:	Q13905
Other Names:	C3G, CRK SH3-binding GNRP, GRF2, RPGF1, Rap guanine nucleotide exchange factor 1, guanine nucleotide-releasing factor 2
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 RapGEF1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RapGEF1 antibody detects endogenous levels of total RapGEF1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
References:	Tanaka S., Proc. Natl. Acad. Sci. U.S.A. 91:3443-3447(1994). Knudsen B., J. Biol. Chem. 269:32781-32787(1994). Humphray S.J., Nature 429:369-374(2004).