

ARHGEF12 Antibody [#C18388]

Catalog Number:	C18388
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NZN5
Other Names:	ARHGC; ARHGEF12; Leukemia-associated RhoGEF; RHO guanine nucleotide exchange factor (GEF) 12 (LARG); Rho guanine nucleotide exchange factor 12
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 ARHGEF12.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ARHGEF12 antibody detects endogenous levels of total ARHGEF12 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Kourlas P.J., Proc. Natl. Acad. Sci. U.S.A. 97:2145-2150(2000). The MGC Project Team, Genome Res. 14:2121-2127(2004). Nagase T., DNA Res. 4:141-150(1997).