

GIMAP2 Antibody [#C16003]

Catalog Number:	C16003
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UG22
Other Names:	gima2; gimap2; gtpase imap family member 2; gtpase, imap family member 2; himap2; imap2; immunity-associated protein 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 GIMAP2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GIMAP2 antibody detects endogenous levels of total GIMAP2 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	The German cDNA consortium; Submitted (AUG-1999) to the EMBL/GenBank/DDBJ databases. The MGC Project Team; Genome Res. 14:2121-2127(2004).