

GIMAP4 Antibody [#C16004]

Catalog Number:	C16004
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NUV9
Other Names:	GIMA4; GTPase, IMAP 4; IAN1; IMAP family member 4; IMAP4; immune associated nucleotide 1; immunity associated protein 4; immunity-associated 4; immunity-associated nucleotide 1 protein; MSTP062
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 GIMAP4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GIMAP4 antibody detects endogenous levels of total GIMAP4 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Cambot M., Blood 99:3293-3301(2002). Liu Y.Q., Submitted (DEC-1998) to the EMBL/GenBank/DDBJ databases. Ota T., Nat. Genet. 36:40-45(2004).