

GNG5 Antibody [#C16046]

Catalog Number:	C16046
Concentration:	1mg/ml
Swiss-Prot No.:	P63218
Other Names:	Guanine nucleotide-binding protein G(I)/G(S)/G(O) subunit gamma-5
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 GNG5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GNG5 antibody detects endogenous levels of total GNG5 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:40000
References:	Liu B., Biochem. Biophys. Res. Commun. 251:88-94(1998). Mao M., Proc. Natl. Acad. Sci. U.S.A. 95:8175-8180(1998). Puhl H.L. III, Submitted (MAR-2002) to the EMBL/GenBank/DBJ databases.