

## GNAT1 Antibody [#C16062]

<b>Catalog Number:</b>	C16062
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P11488
<b>Other Names:</b>	GBT1; GNATR; guanine nucleotide-binding protein G; transducin alpha-1 ch
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human GNAT1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	GNAT1 antibody detects endogenous levels of total GNAT1 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Fong S.-L.; Nucleic Acids Res. 20:2865-2870(1992). Lerea C.L., Neuron 3:367-376(1989). van Dop C., Nucleic Acids Res. 17:4887-4887(1989).