

GABA-RB Antibody [#B0011]

Catalog Number:	B0011
Concentration:	1mg/ml
Swiss-Prot No.:	P18505
Other Names:	GABA(A) receptor, GABA(A)R beta-1, GBRB1, Gamma-aminobutyric-acid receptor beta-1 subunit precursor
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 GABA-RB.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GABA-RB antibody detects endogenous levels of total GABA-RB protein.
Reactivity:	Human, Mouse, Rat
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Schofield P.R., FEBS Lett. 244:361-364(1989). Kirkness E.F., Genomics 10:985-995(1991). Ota T., Nat. Genet. 36:40-45(2004).