

ADRA1B Antibody [#G024]

Catalog Number:	G024
Concentration:	1mg/ml
Swiss-Prot No.:	P35368
Other Names:	Alpha-1B adrenergic receptor; Alpha 1B-adrenoceptor; Alpha 1B-adrenoreceptor; ADRA1B
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 ADRA1B.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ADRA1B antibody detects endogenous levels of total ADRA1B protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 93%, Positives = 93%); Rat (Identities = 93%, Positives = 93%)
Applications:	IF: 1:500~1:1000 ELISA: 1:20000
References:	Ramarao C.S., J. Biol. Chem. 267:21936-21945(1992). Forray C., Mol. Pharmacol. 45:703-708(1994). Schwinn D.A., J. Pharmacol. Exp. Ther. 272:134-142(1995).