

ADRA1A Antibody [#G025]

Catalog Number:	G025
Concentration:	1mg/ml
Swiss-Prot No.:	P35348
Other Names:	Alpha-1A adrenergic receptor; Alpha 1A-adrenoceptor; Alpha 1A-adrenoreceptor; Alpha-1C adrenergic receptor; Alpha-adrenergic receptor 1c; ADRA1A; ADRA1C
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 ADRA1A.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ADRA1A antibody detects endogenous levels of total ADRA1A protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 90%, Positives = 90%); Rat (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:10000
References:	Hirasawa A., Biochem. Biophys. Res. Commun. 195:902-909(1993). Weinberg D.H., Biochem. Biophys. Res. Commun. 201:1296-1304(1994). Forray C., Mol. Pharmacol. 45:703-708(1994).