

5-HT-1B Antibody [#C12009]

Catalog Number:	C12009
Concentration:	1mg/ml
Swiss-Prot No.:	P28222
Other Names:	5-hydroxytryptamine receptor 1B, 5-HT-1B, Serotonin receptor 1B, 5-HT1B, 5-HT-1D-beta, Serotonin 1D beta receptor, S12, HTR1B, HTR1DB
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 5-HT-1B.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	5-HT-1B antibody detects endogenous levels of total 5-HT-1B protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Weinshank R.L., Proc. Natl. Acad. Sci. U.S.A. 89:3630-3634(1992). Demchyshyn L., Proc. Natl. Acad. Sci. U.S.A. 89:5522-5526(1992). Mochizuki D., Biochem. Biophys. Res. Commun. 185:517-523(1992).