

HTR4 Antibody [#G014]

Catalog Number:	G014
Concentration:	1mg/ml
Swiss-Prot No.:	Q13639
Other Names:	5-hydroxytryptamine receptor 4; 5-HT-4; 5-HT4; Serotonin receptor 4; HTR4
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 HTR4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HTR4 antibody detects endogenous levels of total HTR4 protein.
Reactivity:	Human, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
References:	Blondel O., J. Neurochem. 70:2252-2261(1998). Claeyens S., NeuroReport 8:3189-3195(1997). Claeyens S., Mol. Pharmacol. 55:910-920(1999).