

## 5-HT-1A Antibody [#C12008]

<b>Catalog Number:</b>	C12008
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P08908
<b>Other Names:</b>	5-HT1A, 5-hydroxytryptamine 1A receptor, 5HT1A, G-21, HTR1A, Serotonin receptor 1A
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human 5-HT-1A.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	5-HT-1A antibody detects endogenous levels of total 5-HT-1A protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Kobilka B.K., Nature 329:75-79(1987). The MGC Project Team, Genome Res. 14:2121-2127(2004). Parks C.L., Nucleic Acids Res. 19:7155-7160(1991).