

5-HT-3A Antibody [#C12016]

| | |
|---------------------------|--|
| Catalog Number: | C12016 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | P46098 |
| Other Names: | 5-HT-3, 5-HT3R, 5-hydroxytryptamine 3 receptor, 5HT3R, HTR3, HTR3A, serotonin-gated ion channel receptor |
| 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-3A. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | 5-HT-3A antibody detects endogenous levels of total 5-HT-3A protein. |
| Reactivity: | Human |
| Applications: | WB: 1:500~1:1000 ELISA: 1:10000 |
| References: | Miyake A., Mol. Pharmacol. 48:407-416(1995). Belelli D., Mol. Pharmacol. 48:1054-1062(1995). Yamada K., Biol. Psychiatry 60:192-201(2006). |