

OR5A2 Antibody [#G622]

| | |
|---------------------------|--|
| Catalog Number: | G622 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | Q8NGI9 |
| Other Names: | Olfactory receptor 5A2 |
| 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 C-terminal of human OR5A2. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | OR5A2 antibody detects endogenous levels of total OR5A2 protein. |
| Reactivity: | Human |
| Applications: | WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000 |
| References: | Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases. Fuchs T., Genomics 80:295-302(2002). |