

Cytochrome P450 3A4/5 Antibody [#C12276]

| | |
|---------------------------|---|
| Catalog Number: | C12276 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | P05184/P08684/P20815 |
| Other Names: | Cytochrome P450 3A4, EC 1.14.13.67, Quinine 3-monooxygenase, CYP11A4, Nifedipine oxidase, Nifedipine oxidase, Cytochrome P450 3A3, CYP11A3, HLp, Taurochenodeoxycholate 6-alpha-hydroxylase, EC 1.14.13.97, NF-25, P450-PCN1, CYP3A3, CP3A4 |
| 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 Cytochrome P450 3A4/5. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | Cytochrome P450 3A4/5 antibody detects endogenous levels of total Cytochrome P450 3A4/5 protein. |
| Reactivity: | Human |
| Applications: | WB: 1:500~1:1000 ELISA: 1:10000 |
| References: | Chen Q., ghua Yu Fang Yi Xue Za Zhi 32:281-284(1998). Gellner K., Pharmacogenetics 11:111-121(2001). Hsieh K.-P., Drug Metab. Dispos. 29:268-273(2001). |