

## COX5B Antibody [#C12238]

|                           |   |
|---------------------------|---|
| <b>Catalog Number:</b>    | C12238  |
| <b>Concentration:</b>     | 1mg/ml  |
| <b>Swiss-Prot No.:</b>    | P10606  |
| <b>Other Names:</b>       | Cytochrome c oxidase polypeptide Vb, Cytochrome c oxidase subunit 5B, mitochondrial precursor, EC 1.9.3.1   |
| <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 COX5B.  |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.                                   |
| <b>Specificity:</b>       | COX5B antibody detects endogenous levels of total COX5B protein.  |
| <b>Reactivity:</b>        | Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)                     |
| <b>Applications:</b>      | IHC: 1:50~1:100 ELISA: 1:20000  |
| <b>References:</b>        | Lomax M.I., Genomics 10:1-9(1991).<br>Bachman N.J., Arch. Biochem. Biophys. 333:152-162(1996).<br>The MGC Project Team, Genome Res. 14:2121-2127(2004). |