

Cytochrome P450 4X1 Antibody [#C12290]

Catalog Number:	C12290
Concentration:	1mg/ml
Swiss-Prot No.:	Q8N118
Other Names:	CP4X1; CYP1VX1; Cytochrome P450 4X1; EC 1.14.14.1
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 4X1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Cytochrome P450 4X1 antibody detects endogenous levels of total Cytochrome P450 4X1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Clark H.F., Genome Res. 13:2265-2270(2003). Ota T., Nat. Genet. 36:40-45(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004).