

Cytochrome P450 39A1 Antibody [#C12275]

Catalog Number:	C12275
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NYL5
Other Names:	Cytochrome P450 39A1, EC 1.14.13.99, 24-hydroxycholesterol 7-alpha-hydroxylase, Oxysterol 7-alpha-hydroxylase, Hcyp39a1, CP39A
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 39A1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Cytochrome P450 39A1 antibody detects endogenous levels of total Cytochrome P450 39A1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
References:	Li-Hawkins J., . Biol. Chem. 275:16543-16549(2000). The MGC Project Team, ome Res. 14:2121-2127(2004).