

Cytochrome P450 1A1/2 Antibody [#C12249]

Catalog Number:	C12249
Concentration:	1mg/ml
Swiss-Prot No.:	P04798/P05177
Other Names:	AHH, AHRR, CP11, CYP1, CYP1A1, EC 1.14.14.1, P1-450, P450 form 6, P450-C, P450-P1, P450DX, aryl hydrocarbon hydroxylase, cytochrome P1-450, dioxin-inducible, cytochrome P450 1A1, cytochrome P450 1A1 variant, cytochrome P450, family 1, subfamily A, polypeptide 1, cytochrome P450, subfamily I (aromatic compound-inducible), polypeptide 1, flavoprotein-linked monooxygenase
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 1A1/2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Cytochrome P450 1A1/2 antibody detects endogenous levels of total Cytochrome P450 1A1/2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Jaiswal A.K., Nucleic Acids Res. 13:4503-4520(1985). The MGC Project Team, Genome Res. 14:2121-2127(2004). Olsen J.V., II 127:635-648(2006).