

Cytochrome P450 2D6 Antibody [#C21042]

Catalog Number:	C21042
Concentration:	1mg/ml
Swiss-Prot No.:	Q2XND8/P10635
Other Names:	CP2D6; CPD6; CYP2D; CYP2D@; CYP2DL1; CYP1ID6; cytochrome P450 2D6; cytochrome P450, family 2, subfamily D, polypeptide 6; debrisoquine 4-hydroxylase; EC 1.14.14.1; P450-DB1
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 2D6.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Cytochrome P450 2D6 antibody detects endogenous levels of total Cytochrome P450 2D6 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Gonzalez F.J., Genomics 2:174-179(1988). Gonzalez F.J., Nature 331:442-446(1988). Kimura S., Am. J. Hum. Genet. 45:889-904(1989).