

Cytochrome P450 27A1 Antibody [#C12257]

Catalog Number:	C12257
Concentration:	1mg/ml
Swiss-Prot No.:	Q02318
Other Names:	25-hydroxylase, 5-beta-cholestane-3-alpha,7-alpha,12-alpha-triol 27-hydroxylase, CP27A, Cytochrome P-450C27/25, Cytochrome P450 27, mitochondrial, EC 1.14.13.15, Sterol 26-hydroxylase, Sterol 27-hydroxylase, Vitamin D
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 27A1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Cytochrome P450 27A1 antibody detects endogenous levels of total Cytochrome P450 27A1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Cali J.J., J. Biol. Chem. 266:7774-7778(1991). Guo Y.-D., Proc. Natl. Acad. Sci. U.S.A. 90:8668-8672(1993). Leitersdorf E., J. Clin. Invest. 91:2488-2496(1993).