

ATP5H Antibody [#C14591]

Catalog Number:	C14591
Concentration:	1mg/ml
Swiss-Prot No.:	O75947
Other Names:	ATP synthase D chain, mitochondrial; ATP synthase, H ⁺ transporting, mitochondrial F0 complex, subunit d; ATP synthase, H ⁺ transporting, mitochondrial F1F0, subunit d; ATP5JD; ATPQ; EC 3.6.3.14; My032 protein
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 ATP5H.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ATP5H antibody detects endogenous levels of total ATP5H protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Lee H.C., Submitted (AUG-1998) to the EMBL/GenBank/DDBJ databases. Zhang Q.-H., Genome Res. 10:1546-1560(2000). Mao Y.M., Submitted (APR-1998) to the EMBL/GenBank/DDBJ databases.