

ATP5D Antibody [#C14592]

Catalog Number:	C14592
Concentration:	1mg/ml
Swiss-Prot No.:	P30049
Other Names:	ATP synthase subunit delta, mitochondrial
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 ATP5D.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ATP5D antibody detects endogenous levels of total ATP5D protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 83%, Positives = 91%); Rat (Identities = 91%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Grimwood J., Nature 428:529-535(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004). Hochstrasser D.F., Electrophoresis 13:992-1001(1992).