

ATP5G2 Antibody [#C14598]

Catalog Number:	C14598
Concentration:	1mg/ml
Swiss-Prot No.:	Q06055
Other Names:	ATP synthase lipid-binding protein, mitochondrial; ATP synthase proteolipid P2; ATPase protein 9; ATPase subunit c
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 ATP5G2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ATP5G2 antibody detects endogenous levels of total ATP5G2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Dyer M.R., Biochem. J. 293:51-64(1993). Higuti T., Biochim. Biophys. Acta 1173:87-90(1993). Farrell L.B., Biochem. Biophys. Res. Commun. 144:1257-1264(1987).