

COX41 Antibody [#C12235]

Catalog Number:	C12235
Concentration:	1mg/ml
Swiss-Prot No.:	P13073
Other Names:	COX IV-1, COX4, COXIV, EC 1.9.3.1, cytochrome c oxidase polypeptide IV
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 N-terminal of human COX41.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	COX41 antibody detects endogenous levels of total COX41 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Lomax M.I., Gene 86:209-216(1990). Bachman N.J., Mamm. Genome 10:506-512(1999). van Kuilenburg A.B.P., chim. Biophys. Acta 1119:218-224(1992).