

COX7S/A2 Antibody [#C12243]

Catalog Number:	C12243
Concentration:	1mg/ml
Swiss-Prot No.:	O60397
Other Names:	Cytochrome c oxidase subunit VIIa 3, mitochondrial [Precursor], COX7AP2, COX7A3, COX7AL2, COX7S
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 COX7S/A2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	COX7S/A2 antibody detects endogenous levels of total COX7S/A2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	Hillier L.W., Nature 434:724-731(2005).