

## NDUFA4 Antibody [#C16821]

<b>Catalog Number:</b>	C16821
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O00483
<b>Other Names:</b>	CI-MLRQ; complex I-MLRQ; EC 1.6.5.3; EC 1.6.99.3; MLRQ; NADH dehydrogenase (ubiquinone) 1 alpha subcomplex, 4 (9kD, MLRQ); NADH dehydrogenase (ubiquinone) 1 alpha subcomplex, 4, 9kDa; NADH dehydrogenase [ubiquinone] 1 alpha subcomplex 4; NADH-ubiquinone oxidoreductase MLRQ; NDUA4
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human NDUFA4.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	NDUFA4 antibody detects endogenous levels of total NDUFA4 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 92%); Rat (Identities = 92%, Positives = 92%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Kim J.W., Biochem. Mol. Biol. Int. 43:669-675(1997). Kanagarajah D., Submitted (NOV-1999) to the EMBL/GenBank/DDBJ databases. Ebert L., Submitted (MAY-2004) to the EMBL/GenBank/DDBJ databases.