

NDUC2 Antibody [#C10222]

Catalog Number:	C10222
Concentration:	1mg/ml
Swiss-Prot No.:	O95298
Other Names:	NADH dehydrogenase [ubiquinone] 1 subunit C2, EC 1.6.5.3, EC 1.6.99.3, NADH-ubiquinone oxidoreductase subunit B14.5b, Complex I-B14.5b, CI-B14.5b, Lung cancer oncogene 1 protein, HLC-1, NDUFC2, HLC1
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 NDUC2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NDUC2 antibody detects endogenous levels of total NDUC2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	Loeffen J.L.C.M., Biochem. Biophys. Res. Commun. 253:415-422(1998). Zhang Q.-H., Genome Res. 10:1546-1560(2000). Wiemann S., Genome Res. 11:422-435(2001).