

## NDUFB10 Antibody [#C16829]

<b>Catalog Number:</b>	C16829
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O96000
<b>Other Names:</b>	EC 1.6.5.3; EC 1.6.99.3; NADH-ubiquinone oxidoreductase PDSW; NDUBA
<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 NDUFB10.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	NDUFB10 antibody detects endogenous levels of total NDUFB10 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Loeffen J.L.C.M., Biochem. Biophys. Res. Commun. 253:415-422(1998). Wang L., Submitted (AUG-1998) to the EMBL/GenBank/DDBJ databases. Zhang Q.-H., Genome Res. 10:1546-1560(2000).