

NDUFV2 Antibody [#C16842]

Catalog Number:	C16842
Concentration:	1mg/ml
Swiss-Prot No.:	P19404
Other Names:	EC 1.6.5.3; EC 1.6.99.3; NDUV2
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 NDUFV2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NDUFV2 antibody detects endogenous levels of total NDUFV2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Pilkington S.J., Biochemistry 28:3257-3264(1989). The MGC Project Team; Genome Res. 14:2121-2127(2004).