

PNPT1 Antibody [#C17774]

Catalog Number:	C17774
Concentration:	1mg/ml
Swiss-Prot No.:	Q8TCS8
Other Names:	Polyribonucleotide nucleotidyltransferase 1, mitochondrial; EC 2.7.7.8; PNPase 1; Polynucleotide phosphorylase-like protein; PNPase old-35; 3'-5' RNA exonuclease OLD35;3'-5' RNA exonuclease OLD35; EC 2.7.7.8; PNPASE; PNPase 1; PNPase old-35; PNPT1; polynucleotide phosphorylase-like protein; polyribonucleotide nucleotidyltransferase 1, mitochondrial
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 PNPT1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PNPT1 antibody detects endogenous levels of total PNPT1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Raijmakers R.,J. Mol. Biol. 323:653-663(2002). Leszczyniecka M.,Proc. Natl. Acad. Sci. U.S.A. 99:16636-16641(2002). Hillier L.W., Nature 434:724-731(2005).