

PINX1 Antibody [#C11479]

Catalog Number:	C11479
Concentration:	1mg/ml
Swiss-Prot No.:	Q96BK5
Other Names:	Pin2-interacting protein X1, TRF1-interacting protein 1, Liver-related putative tumor suppressor, Protein 67-11-3, PINX1, LPTL, LPTS
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 PINX1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PINX1 antibody detects endogenous levels of total PINX1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Liao C., Hepatology 32:721-727(2000). Zhou X.Z., Cell 107:347-359(2001). Song H., Biochem. Biophys. Res. Commun. 316:1116-1123(2004).