

CLPT1 Antibody [#G070]

Catalog Number:	G070
Concentration:	1mg/ml
Swiss-Prot No.:	O96005
Other Names:	Cleft lip and palate transmembrane protein 1; CLPTM1
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 CLPT1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CLPT1 antibody detects endogenous levels of total CLPT1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Yoshiura K., Genomics 54:231-240(1998). Totoki Y., Submitted (APR-2005) to the EMBL/GenBank/DDBJ databases. Lewandrowski U., Mol. Cell. Proteomics 5:226-233(2006).