

GLB1L3 Antibody [#C14733]

Catalog Number:	C14733
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NCI6
Other Names:	Beta-galactosidase-1-like protein 3; EC 3.2.1.-
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 GLB1L3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GLB1L3 antibody detects endogenous levels of total GLB1L3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Ota T., Nat. Genet. 36:40-45(2004). Taylor T.D., Nature 440:497-500(2006). The MGC Project Team; Genome Res. 14:2121-2127(2004).