

GPR137C Antibody [#G264]

Catalog Number:	G264
Concentration:	1mg/ml
Swiss-Prot No.:	Q8N3F9
Other Names:	Integral membrane protein GPR137C; Transmembrane 7 superfamily member 1-like 2 protein; GPR137C; TM7SF1L2
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 GPR137C.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR137C antibody detects endogenous levels of total GPR137C protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Heilig R., Nature 421:601-607(2003). Vassilatis D.K., Proc. Natl. Acad. Sci. U.S.A. 100:4903-4908(2003). The German cDNA consortium; Submitted (SEP-2004) to the EMBL/GenBank/DDBJ databases.