

ISL2 Antibody [#C11470]

Catalog Number:	C11470
Concentration:	1mg/ml
Swiss-Prot No.:	Q96A47
Other Names:	Insulin gene enhancer protein ISL-2, Islet-2, ISL2
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 ISL2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ISL2 antibody detects endogenous levels of total ISL2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	The MGC Project Team, Genome Res. 14:2121-2127(2004). Beausoleil S.A., Proc. Natl. Acad. Sci. U.S.A. 101:12130-12135(2004).