

INSL4 Antibody [#C15598]

Catalog Number:	C15598
Concentration:	1mg/ml
Swiss-Prot No.:	Q14641
Other Names:	Early placenta insulin-like peptide [Precursor]; EPIL; Placentin; Insulin-like peptide 4; Early placenta insulin-like peptide B chain; Early placenta insulin-like peptide A chain; INSL4
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 INSL4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	INSL4 antibody detects endogenous levels of total INSL4 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Chassin D., Genomics 29:465-470(1995). Humphray S.J., Nature 429:369-374(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004).