

CIDEC Antibody [#C12188]

Catalog Number:	C12188
Concentration:	1mg/ml
Swiss-Prot No.:	Q96AQ7
Other Names:	Cell death activator CIDE-3, Cell death-inducing DFFA-like effector protein C, Fat-specific protein FSP27 homolog, FSP27
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 CIDEC.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CIDEC antibody detects endogenous levels of total CIDEC protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Liang L., Biochem. J. 370:195-203(2003). Ota T., Nat. Genet. 36:40-45(2004). The MGC Project Team, Genome Res. 14:2121-2127(2004).