

DNL4 Antibody [#C13046]

Catalog Number:	C13046
Concentration:	1mg/ml
Swiss-Prot No.:	P49917
Other Names:	DNA ligase 4, EC 6.5.1.1, DNA ligase IV, Polydeoxyribonucleotide synthase [ATP] 4, LIG4
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 DNL4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	DNL4 antibody detects endogenous levels of total DNL4 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Wei Y.-F., Mol. Cell. Biol. 15:3206-3216(1995). Dunham A., Nature 428:522-528(2004). Robins P., J. Biol. Chem. 271:24257-24261(1996).