

GNL3L Antibody [#C16072]

Catalog Number:	C16072
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NVN8
Other Names:	guanine nucleotide binding protein-like 3 (nucleolar)-like; guanine nucleotide-binding protein-like 3-like protein
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 GNL3L.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GNL3L antibody detects endogenous levels of total GNL3L protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Ota T., Nat. Genet. 36:40-45(2004). Ross M.T., Nature 434:325-337(2005). The MGC Project Team; Genome Res. 14-2121-2127(2004).