

GNL1 Antibody [#C16070]

Catalog Number:	C16070
Concentration:	1mg/ml
Swiss-Prot No.:	P36915
Other Names:	GTP-binding protein HSR1; guanine nucleotide binding protein-like 1; guanine nucleotide-binding protein-like 1; HSR1; HSR1 GTP-binding 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 GNL1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GNL1 antibody detects endogenous levels of total GNL1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Vernet C., Mamm. Genome 5:100-105(1994). Kalnine N., Submitted (MAY-2003) to the EMBL/GenBank/DDBJ databases. Mungall A.J., Nature 425:805-811(2003).