

NSG1 Antibody [#C17009]

Catalog Number:	C17009
Concentration:	1mg/ml
Swiss-Prot No.:	P42857
Other Names:	brain neuron cytoplasmic protein 1; d4s234e; neuron-specific protein family member 1; NSG1
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 NSG1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NSG1 antibody detects endogenous levels of total NSG1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Carlock L., Brain Res. Mol. Brain Res. 42:202-212(1996). The MGC Project Team; Genome Res. 14:2121-2127(2004).