

NSG2 Antibody [#C17010]

Catalog Number:	C17010
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y328
Other Names:	Neuron-specific protein family member 2, Protein p19; HMP19
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 NSG2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NSG2 antibody detects endogenous levels of total NSG2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Jin J., Proc. Natl. Acad. Sci. U.S.A. 97:9543-9548(2000). The MGC Project Team; Genome Res. 14:2121-2127(2004).