

GSC2 Antibody [#C10085]

Catalog Number:	C10085
Concentration:	1mg/ml
Swiss-Prot No.:	O15499
Other Names:	Homeobox protein goosecoid-2, GSC-2, Homeobox protein goosecoid-like, GSCL
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 GSC2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GSC2 antibody detects endogenous levels of total GSC2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Funke B., Genomics 46:364-372(1997). Gottlieb S., Am. J. Hum. Genet. 60:1194-1201(1997).