

CSPG5 Antibody [#C15097]

Catalog Number:	C15097
Concentration:	1mg/ml
Swiss-Prot No.:	O95196
Other Names:	Chondroitin sulfate proteoglycan 5; Neuroglycan C; Acidic leucine-rich EGF-like domain-containing brain 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 CSPG5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CSPG5 antibody detects endogenous levels of total CSPG5 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Yasuda Y., Neurosci. Res. 32:313-322(1998). Aono S., Submitted (DEC-2001) to the EMBL/GenBank/DDBJ databases. Muzny D.M., Nature 440:1194-1198(2006).