

CSGALNACT1 Antibody [#C15093]

Catalog Number:	C15093
Concentration:	1mg/ml
Swiss-Prot No.:	Q8TDX6
Other Names:	Chondroitin beta-1,4-N-acetylgalactosaminyltransferase 1; EC 2.4.1.174; Beta4GalNAcT-1
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 CSGALNACT1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CSGALNACT1 antibody detects endogenous levels of total CSGALNACT1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%); Rat (Identities = 92%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Uyama T., J. Biol. Chem. 277:8841-8846(2002). Gotoh M., J. Biol. Chem. 277:38189-38196(2002). Clark H.F., Genome Res. 13:2265-2270(2003).