

CHST6 Antibody [#C14929]

Catalog Number:	C14929
Concentration:	1mg/ml
Swiss-Prot No.:	Q9GZX3
Other Names:	Carbohydrate sulfotransferase 6; EC 2.8.2.-; N-acetylglucosamine 6-O-sulfotransferase 5; GlcNAc6ST-5; Corneal N-acetylglucosamine-6-O-sulfotransferase; C-GlcNAc6ST; hCGn6ST; Galactose/N-acetylglucosamine/N-acetylglucosamine 6-O-sulfotransferase 4-beta; GST4-beta
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 CHST6.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CHST6 antibody detects endogenous levels of total CHST6 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:40000
References:	Akama T.O., Nat. Genet. 26:237-241(2000). Hemmerich S., Glycobiology 11:75-87(2001). Bartes A., Biochem. Biophys. Res. Commun. 282:928-933(2001).