

CHST13 Antibody [#C14924]

Catalog Number:	C14924
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NET6
Other Names:	Carbohydrate sulfotransferase 13; EC 2.8.2.5; Chondroitin 4-O-sulfotransferase 3; Chondroitin 4-sulfotransferase 3; C4ST3; C4ST-3
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 CHST13.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CHST13 antibody detects endogenous levels of total CHST13 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Kang H.-G., J. Biol. Chem. 277:34766-34772(2002).