

B3GALT2 Antibody [#C14713]

Catalog Number:	C14713
Concentration:	1mg/ml
Swiss-Prot No.:	O43825
Other Names:	Beta-1,3-galactosyltransferase 2; EC 2.4.1.-; Beta-1,3-GalTase 2; Beta3Gal-T2; UDP-galactose:2-acetamido-2-deoxy-D-glucose 3beta-galactosyltransferase 2
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 B3GALT2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	B3GALT2 antibody detects endogenous levels of total B3GALT2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:60000
References:	Sood R., Genomics 73:211-222(2001). Gregory S.G., Nature 441:315-321(2006). Amado M., J. Biol. Chem. 273:12770-12778(1998).