

POFUT1 Antibody [#C15897]

Catalog Number:	C15897
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H488
Other Names:	EC 2.4.1.221; GDP-fucose protein O-fucosyltransferase 1; KIAA0180; O-FucT-1; OFUT1; peptide-O-fucosyltransferase
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 POFUT1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	POFUT1 antibody detects endogenous levels of total POFUT1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	Wang Y., J. Biol. Chem. 276:40338-40345(2001). Nagase T., DNA Res. 3:17-24(1996). Nakajima D., DNA Res. 9:99-106(2002).