

LFNG Antibody [#C14718]

Catalog Number:	C14718
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NES3
Other Names:	beta-1,3-N-acetylglucosaminyltransferase lunatic fringe; EC 2.4.1.222; LFNG O-fucosylpeptide 3-beta-N-acetylglucosaminyltransferase; O-fucosylpeptide 3-beta-N-acetylglucosaminyltransferase; SCDO3
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 LFNG.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LFNG antibody detects endogenous levels of total LFNG protein.
Reactivity:	Human (Identities = 100%, Positives = 100%) Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 92%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Johnston S.H., Development 124:2245-2254(1997). Cole S.E., Dev. Cell 3:75-84(2002). Shimizu K., J. Biol. Chem. 276:25753-25758(2001).