

GHRHR Antibody [#G271]

Catalog Number:	G271
Concentration:	1mg/ml
Swiss-Prot No.:	Q02643
Other Names:	GHRFR; GHRH receptor; GHRHRpsv; GRF receptor; GRFR; growth hormone-releasing hormone receptor
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 GHRHR.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GHRHR antibody detects endogenous levels of total GHRHR protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Mayo K.E.; Mol. Endocrinol. 6:1734-1744(1992). Gaylinn B.D., Mol. Endocrinol. 7:77-84(1993). Hillier L.W., Nature 424:157-164(2003).