

## LSHR Antibody [#G380]

<b>Catalog Number:</b>	G380
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P22888
<b>Other Names:</b>	Lutropin-choriogonadotropic hormone receptor; LH/CG-R; LSH-R; Luteinizing hormone receptor; LHR; LHCGR; LCGR; LGR2; LHRHR
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human LSHR.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	LSHR antibody detects endogenous levels of total LSHR protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 86%, Positives = 93%); Rat (Identities = 93%, Positives = 100%)
<b>Applications:</b>	IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Minegish T., Biochem. Biophys. Res. Commun. 172:1049-1054(1990). Jia X.-C., Mol. Endocrinol. 5:759-768(1991). Frazier A.L., Mol. Endocrinol. 4:1264-1276(1990).