

## GPR107 Antibody [#G279]

<b>Catalog Number:</b>	G279
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q5VW38
<b>Other Names:</b>	GP107; Lung seven transmembrane receptor 1; Protein GPR107
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 GPR107.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	GPR107 antibody detects endogenous levels of total GPR107 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:5000
<b>References:</b>	Edgar A.J., J. Anat. 200:202-202(2002). Ota T., Nat. Genet. 36:40-45(2004). Humphray S.J., Nature 429:369-374(2004).