

## GPR18 Antibody [#G326]

<b>Catalog Number:</b>	G326
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q14330
<b>Other Names:</b>	N-arachidonyl glycine receptor; NAGly receptor; G-protein coupled receptor 18; GPR18; GPCRW
<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 GPR18.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	GPR18 antibody detects endogenous levels of total GPR18 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Gantz I., Genomics 42:462-466(1997). Kohno M., Biochem. Biophys. Res. Commun. 347:827-832(2006). Dunham A., Nature 428:522-528(2004).