

## FPRL2 Antibody [#G252]

<b>Catalog Number:</b>	G252
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P25089
<b>Other Names:</b>	FMLP-related receptor II;FMLP-R-II; Formylpeptide receptor-like 2; FPRL2;FPRH1
<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 FPRL2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	FPRL2 antibody detects endogenous levels of total FPRL2 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Bao L., Genomics 13:437-440(1992). Durstin M., Biochem. Biophys. Res. Commun. 201:174-179(1994). Kopatz S.A., Submitted (MAR-2003) to the EMBL/GenBank/DDBJ databases.