

FFAR1 Antibody [#G248]

Catalog Number:	G248
Concentration:	1mg/ml
Swiss-Prot No.:	O14842
Other Names:	free fatty acid receptor 1; G-protein coupled receptor 40; GPR40
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 humanFFAR1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	FFAR1antibodydetects endogenous levels of totalFFAR1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000IF: 1:500~1:1000 ELISA: 1:10000
References:	Sawzdargo M., Biochem. Biophys. Res. Commun. 239:543-547(1997). The MGC Project Team; Genome Res. 14:2121-2127(2004). Briscoe C.P., J. Biol. Chem. 278:11303-11311(2003).