

## AGTR1 Antibody [#G212]

<b>Catalog Number:</b>	G212
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P30556
<b>Other Names:</b>	AG2R; AGTR1; AGTR1A; AGTRA; AT1A; AT1AR; AT2R1; Type-1 angiotensin II receptor; Type-1A angiotensin II receptor
<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 AGTR1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	AGTR1 antibody detects endogenous levels of total AGTR1 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Mauzy C.A., Biochem. Biophys. Res. Commun. 186:277-284(1992). Furuta H., Biochem. Biophys. Res. Commun. 183:8-13(1992). Bergsma D.J., Biochem. Biophys. Res. Commun. 183:989-995(1992).