

## PE2R3 Antibody [#G098]

<b>Catalog Number:</b>	G098
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P43115
<b>Other Names:</b>	EP3; EP3-I; EP3-II; EP3-III; EP3-IV; EP3e; PE2R3; PGE receptor, EP3; PGE2-R; prostaglandin E2 receptor EP3
<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 PE2R3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PE2R3 antibody detects endogenous levels of total PE2R3 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: 1:5000
<b>References:</b>	Adam M., FEBS Lett. 338:170-174(1994). Schmid A., Eur. J. Biochem. 228:23-30(1995). Yang J., Biochem. Biophys. Res. Commun. 198:999-1006(1994).