

PE2R4 Antibody [#G099]

Catalog Number:	G099
Concentration:	1mg/ml
Swiss-Prot No.:	P35408
Other Names:	Prostaglandin E2 receptor EP4 subtype; PGE receptor, EP4 subtype; Prostanoid EP4 receptor; PTGER4; PTGER2
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 human PE2R4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PE2R4 antibody detects endogenous levels of total PE2R4 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Bastien L., J. Biol. Chem. 269:11873-11877(1994). An S., Biochem. Biophys. Res. Commun. 197:263-270(1993). Foord S.M., Genomics 35:182-188(1996).