

## OPRK1 Antibody [#G491]

<b>Catalog Number:</b>	G491
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P41145
<b>Other Names:</b>	kappa-type opioid receptor; MSL-1; OPRK; OPRK1; ROR-D
<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 OPRK1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	OPRK1 antibody detects endogenous levels of total OPRK1 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Mansson E., Biochem. Biophys. Res. Commun. 202:1431-1437(1994). Simonin F., Proc. Natl. Acad. Sci. U.S.A. 92:7006-7010(1995). Zhu J., Life Sci. 56:PL201-PL207(1995).