

OPRD1 Antibody [#G490]

Catalog Number:	G490
Concentration:	1mg/ml
Swiss-Prot No.:	P41143
Other Names:	delta-opioid receptor; delta-type opioid receptor; K56; MSL-2; opioid receptor delta; OPRD; OPRD1
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 OPRD1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OPRD1 antibody detects endogenous levels of total OPRD1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Knapp R.J., Life Sci. 54:PL463-PL469(1994). Simonin F., Mol. Pharmacol. 46:1015-1021(1994). Gregory S.G., Nature 441:315-321(2006).