

## PKD1/PKC $\mu$ (Ab-910) Antibody [#B0721]

<b>Catalog Number:</b>	B0721
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q15139
<b>Other Names:</b>	EC 2.7.11.13, KPCD1, PKC-mu, PKCM, PKD, PRKCM, PRKD1, Protein kinase C, mu type, Protein kinase D, kinase PKD1, nPKC-mu
<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 non-phosphopeptide derived from human PKD1/PKC $\mu$ around the phosphorylation site of serine 910 (R-V-S <sup>P</sup> -I-L).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PKD1/PKC $\mu$ (Ab-910) detects endogenous levels of total PKD1/PKC $\mu$ protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Lynn Pritchard, Hum. Mol. Genet., Nov 2000; 9: 2617 - 2627. Klaus B. Piontek, J. Am. Soc. Nephrol., Dec 2004; 15: 3035 - 3043. Saori Nishio, J. Clin. Invest., Apr 2005; 115: 910 - 918. Guanqing Wu, Hum. Mol. Genet., Aug 2002; 11: 1845 - 1854.