

## PKC $\delta$ (Ab-505) Antibody [#B0717]

<b>Catalog Number:</b>	B0717
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q05655
<b>Other Names:</b>	EC 2.7.11.13, KPCD, PKC-delta, PRKCD, kinase PKC-delta, nPKC-delta
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 PKC $\delta$ around the phosphorylation site of threonine 505 (A-S-T <sup>P</sup> -F-C).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PKC $\delta$ (Ab-505) antibody detects endogenous levels of total PKC $\delta$ protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
<b>References:</b>	J. J. Peluso, Endocrinology, Oct 2001; 142: 4203. Peter Hampson, Blood, Aug 2005; 106: 1362 - 1368. K Kitamura, Genes Cells, Apr 2003; 8: 311 - 324.