

## PLCXD1 Antibody [#C17702]

<b>Catalog Number:</b>	C17702
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9NUJ7
<b>Other Names:</b>	Phosphatidylinositol-specific phospholipase C X domain-containing protein 1; PI-PLC X domain-containing protein 1
<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 peptide derived from internal of human PLCXD1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PLCXD1 antibody detects endogenous levels of total PLCXD1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100   ELISA: 1:40000
<b>References:</b>	Ota T., Nat. Genet. 36:40-45(2004). Skaletsky H., Nature 423:825-837(2003). Ross M.T., Nature 434:325-337(2005).