

PKD2 (Ab-876) Antibody [#B0808]

Catalog Number:	B0808
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BZL6
Other Names:	EC 2.7.11.13, HSPC187, KPCD2, KPCO, PRKD2, Protein kinase C, D2 type, Protein kinase D2, Protein kinase D2 homolog, kinase PKD2, nPKC-D2
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 non-phosphopeptide derived from human PKD2 around the phosphorylation site of serine 876 (R-I-S ^P -V-L).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PKD2 (Ab-876) antibody detects endogenous levels of total PKD2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Nicholas Obermüller, Am J Physiol Renal Physiol, Dec 1999; 277: 914. Rong Ma, Mol. Cell. Biol., Sep 2005; 25: 8285 - 8298. Dana R. Rundle, J. Biol. Chem., Jul 2004; 279: 29728 - 29739. Sabine Sturany, J. Biol. Chem., Aug 2002; 277: 29431 - 29436.