

PKC ζ (Ab-560) Antibody [#B0805]

Catalog Number:	B0805
Concentration:	1mg/ml
Swiss-Prot No.:	Q05513
Other Names:	EC 2.7.11.13, KPCZ, PKC-zeta, PKC2, PRKCZ, Protein kinase C zeta type, kinase PKC-zeta, nPKC-zeta
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 PKC ζ around the phosphorylation site of threonine 560 (Q-L-T ^P -P-D).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PKC ζ (Ab-560) antibody detects endogenous levels of total PKC ζ protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Richard Magid, Circ. Res., Sep 2005; 97: 443 - 449. M Fujihara, J. Immunol., Feb 1994; 152: 1898. A. Banan, Am J Physiol Gastrointest Liver Physiol, May 2002; 282: 794. A. Banan, Am J Physiol Gastrointest Liver Physiol, Oct 2002; 283: 909