

PKC α/β (Ab-638) Antibody [#B0716]

Catalog Number:	B0716
Concentration:	1mg/ml
Swiss-Prot No.:	P17252
Other Names:	EC 2.7.11.13, KPCA, PKC III, PKC-A, PKC-III, PKC-alpha, PRKACA, PRKCA, Protein kinase C, alpha type, kinase PKC-alpha
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 α/β around the phosphorylation site of threonine 638 (V-L-T ^P -P-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PKC α/β (Ab-638) antibody detects endogenous levels of total PKC α/β protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Maria Nikodemova, J. Biol. Chem., May 2007; 282: 15208 - 15216. Pedro J. Cejas, J. Biol. Chem., Aug 2005; 280: 28412 - 28423. Dan Chen, J. Biol. Chem., Feb 2004; 279: 4829 - 4839.