

Akt (Ab-450) Antibody [#B0406]

Catalog Number:	B0406
Concentration:	1mg/ml
Swiss-Prot No.:	P31749
Other Names:	AKT, AKT1 kinase, C-AKT, PKB, PKB-alpha, Protein kinase B, RAC, RAC-PK-alpha, RAC-alpha serine/threonine kinase, kinase Akt1
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 Akt around the phosphorylation site of threonine 450 (T-I-T ^P -P-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Akt (Ab-450) antibody detects endogenous levels of total Akt protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Payaningal R. Somanath, J. Biol. Chem., Jun 2007; 10.1074. Xiaoming Ju, PNAS, May 2007; 104: 7438 - 7443. Brian DeBosch, Circulation, May 2006; 113: 2097 - 2104.