

Akt (Ab-326) Antibody [#B0817]

Catalog Number:	B0817
Concentration:	1mg/ml
Swiss-Prot No.:	P31749
Other Names:	AKT, AKT1 kinase, C-AKT, EC 2.7.11.1, 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 tyrosine 326 (N-D-Y ^P -G-R).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Akt (Ab-326) antibody detects endogenous levels of total Akt protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Amy D. Sinor, J. Neurosci., Sep 2004; 24: 8531 - 8541. Daniela Matei, Clin. Cancer Res., Jan 2004; 10: 681. Jesika Faridi, Clin. Cancer Res., Aug 2003; 9: 2933.