

Akt (Ab-129) Antibody [#B0609]

Catalog Number:	B0609
Concentration:	1mg/ml
Swiss-Prot No.:	P31749
Other Names:	AKT; AKT1 kinase; C-AKT; EC 2.7.11.1; kinase Akt1; PKB; PKB-alpha; Protein kinase B; RAC; RAC-alpha serine/threonine kinase; RAC-PK-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 Akt around the phosphorylation site of serine 129 (D-N-S ^P -G-A).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Akt (Ab-129) 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: 1:40000
References:	Jones P.F., Proc. Natl. Acad. Sci. U.S.A. 88:4171-4175(1991). Matsubara A., Diabetologia 44:910-913(2001). The MGC Project Team; Genome Res. 14:2121-2127(2004).