



## Akt(Ab-308) Antibody

Catalog Number 21055

(For Research Use Only)

- Catalog Number:** 21055-1, 21055-2, 21055  
**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl  
**Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg<sup>2+</sup> and Ca<sup>2+</sup>), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.  
**Storage/Stability:** Store at -20°C/1 year  
**Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human Akt around the phosphorylation site of Threonine 308.  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity:** Akt(Ab-308) Antibody detects endogenous levels of total Akt protein.  
**Reactivity:** Human, Mouse, Rat  
**Applications:** WB: 1:500~1:1000 IHC1:50~1: 100 ELISA: 1:4000  
**Swiss-Prot No. :** P31749  
**References:** Tremblay F, et al. (2005)Diabetes; 54(9): 2674-84.  
Xu BE, et al. (2005)J Biol Chem; 280(40): 34218-23.  
Samuels Y, et al. (2005)Cancer Cell; 7(6): 561-73.  
Di Maira G, et al. (2005)Cell Death Differ; 12(6): 668-77.

