

Akt (Ab-308) Antibody [#B7005]

- Catalog Number:** B7005-1, B7005-2, B7005
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), 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 IHC: 1:50~1:100
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.