

Akt(Ab-473) Antibody [#B7004]

- Catalog Number:** B7004-1, B7004-2, B7004
- 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:** This antiserum was produced against synthesized non-phosphopeptide derived from human Akt around the phosphorylation site of Serine 473.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** Akt(Ab-473) Antibody detects endogenous levels of total Akt protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:100~1:500 IHC: 1:50~1:100
ELISA: 1:10000
- Swiss-Prot No. :** P31749
- References:** Baudhuin LM, et al. (2004) FASEB J Feb; 18(2): 341-3.
Min YH, et al. (2004) Cancer Res; 64(15): 5225-31.
Feng J, et al. (2004) J Biol Chem; 279(34): 35510-7.
Ayala G, et al. (2004) Clin Cancer Res; 10(19): 6572-8.
Lungu AO, et al. (2004) J Biol Chem; 279(47): 48794-800.