

Akt2 (Ab-474) Antibody [#B7006]

- Catalog Number:** B7006-1, B7006-2, B7006
- 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 Akt2 around the phosphorylation site of serine 474.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** Akt2 (Ab-474) antibody detects endogenous levels of total Akt2 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:10000
- Swiss-Prot No. :** P31751
- References:** Sun M, et al. (2001) Cancer Res; 61(16): 5985-91.
Yuan ZQ, et al. (2000) Oncogene; 19(19): 2324-30.
Meier R, et al. (1997) J Biol Chem; 272(48): 30491-7.