

Instruction Manual

| | |
|---------------------|---|
| Catalog Number | PK-AB718-3519P |
| Quantity | 50 µg |
| Source | 16 amino acids near the amino terminus of human Akt |
| Formulation | Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH ₂ PO ₄ , 10 mM Na ₂ HPO ₄ , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide. |
| Reconstitution | During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap. |
| Storage & Stability | Store Akt1 peptide at -20°C, stable for one year. |
| Application | Akt1 peptide is used for blocking the activity of the Akt antibody. |
| References | <p>Bellacosa A, Testa JR, Staal SP, et al. A retroviral oncogene, akt, encoding a serine-threonine kinase containing an SH2-like region. <i>Science</i> 1991; 254:274-7.</p> <p>Wagner H. The immunobiology of the TLR9 subfamily. <i>Trends Immunol.</i> 2004; 381-6.</p> <p>Nishiya T and DeFranco AL. Ligand-regulated chimeric receptor approach reveals distinctive subcellular localization and signaling properties of the Toll-like receptors. <i>J. Biol. Chem.</i> 2004; 279:19008-17.</p> <p>Zhang G and Ghosh S. Negative regulation of Toll-like receptor-mediated signaling by Tollip. <i>J. Biol. Chem.</i> 2002; 277:7059-65.</p> |

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.
