

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3779P
<b>Quantity</b>	50 µg
<b>Source</b>	16 amino acids near the amino terminus of human β-Actin
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide.
<b>Reconstitution</b>	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.
<b>Storage &amp; Stability</b>	Store Beta-actin peptide at -20°C, stable for one year.
<b>Application</b>	beta-actin peptide is used for blocking the activity of beta-actin antibody.
<b>References</b>	Lambrechts A, Van Troys, M and Ampe C. The actin cytoskeleton in normal and pathological cell motility. <i>Int. J. Biochem. Cell Biol.</i> 2004; 36:1890-909. Yang RB, Mark MR, Gurney AL, et al. Signaling events induced by lipopolysaccharide-activated toll-like receptor 2. <i>J. Immunol.</i> 1999; 163:639-43. Hardy MP and O'Neill LAJ. The murine IRAK2 gene encodes four alternately spliced isoforms, two of which are inhibitory. <i>J. Biol. Chem.</i> 2004; 279:27699-708. van Eyndhoven WG, Gamper CJ, Cho E, et al. TRAF-3 mRNA splice-deletion variants encode isoforms that induce NF0kappaB activation. <i>Mol. Immunol.</i> 1999; 36:647-58.

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