

Instruction Manual

Catalog Number	PK-AB718-3777P
Quantity	50 µg
Source	16 amino acids near the carboxy terminus of human β-Actin
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 Beta-actin peptide at -20°C, stable for one year.
Application	beta-actin peptide is used for blocking the activity of b-actin antibody.
References	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.

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