

Instruction Manual

Catalog Number	PK-AB718-3885P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of human PAK2
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 PAK2 peptide at -20°C, stable for one year.
Application	PAK2 peptide is used for blocking the activity of PAK2 antibody.
References	Jaffer ZM and Chernoff J. p21-activated kinases: three more join the Pak. <i>Int. J. Biochem. Cell Biol.</i> 2002; 34:713-7. Cato ACB and Mink S. BAG-1 family of cochaperones in the modulation of nuclear receptor action. <i>J. Steroid Biochem. & Mol. Biol.</i> 2001; 78:379-88. Kajewska M, Turner BC, Shabaik A, et al. Expression of BAG-1 protein correlates with aggressive behavior of prostate cancers. <i>Prostate</i> 2006; 66:801-10.

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