

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-4085P
<b>Quantity</b>	50 µg
<b>Source</b>	23 amino acids near the carboxy terminus of human BRSK2
<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 BRSK2 peptide at -20°C, stable for one year.
<b>Application</b>	BRSK2 peptide is used for blocking the activity of BRSK2 antibody.
<b>References</b>	Manning G, Whyte DB, Martinez R, et al. The protein kinase complement of the human genome. Science 2002; 298:1912-34.

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