

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-4299P
<b>Quantity</b>	50 µg
<b>Source</b>	15 amino acids near the amino terminus of human PTK7
<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 PTK7 peptide at -20°C, stable for one year.
<b>Application</b>	PTK7 peptide is used for blocking the activity of PTK7 antibody.
<b>References</b>	Park SK, Lee HS and Lee ST. Characterization of the human full-length PTK7 cDNA encoding a receptor protein kinase-like molecule closely related to chick KLG. J. Biochem. 1996; 119:235-9.

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