

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3321P
<b>Quantity</b>	50 µg
<b>Source</b>	15 amino acids near the carboxy terminus of human PIKE
<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 PIKE peptide at -20°C, stable for one year.
<b>Application</b>	PIKE peptide is used for blocking the activity of Anit-PIKE antibody.
<b>References</b>	<p>Ye K, Hurt KJ, Wu FY, et al. Pike. A nuclear gtpase that enhances PI3kinase activity and is regulated by protein 4.1N. Cell 2000; 103:919-30.</p> <p>Hatzioannou T, Perez-Caballero D, Yang A, et al. Retrovirus resistance factors REF1 and Lv1 are species-specific variants of TRIM5alpha. Proc. Nat'l. Acad. Sci. USA 2004; 101:10774-9</p> <p>Xu L, Yang L, Moitra PK, et al. BTBD1 and BTBD2 colocalize to cytoplasmic bodies with the RBCC/tripartite motif protein, TRIM5 . Exp. Cell Res. 2003; 288:84-93.</p>

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