

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-4419P
<b>Quantity</b>	50 µg
<b>Source</b>	17 amino acids near the carboxy terminus of human Plxdc2
<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 Plxdc2 peptide at -20°C, stable for one year.
<b>Application</b>	Plxdc2 peptide is used for blocking the activity of Plxdc2 antibody.
<b>References</b>	Carson-Weber EB, Watkins DN, Nanda A, et al. Cell surface tumor epithelial markers are conserved in mice and humans. Cancer Res. 2001; 61:6649-55.

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