

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3151P
<b>Quantity</b>	50 µg
<b>Source</b>	Amino acids near the carboxy terminus of human PHAP I
<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 PHAP I peptide at -20°C, stable for one year.
<b>Application</b>	PHAP I peptide is used for blocking the activity of PHAP I antibody.
<b>References</b>	Jiang X, Kim HE, Shu H, Zhao Y, Zhang H, Kofron J, Donnelly J, Burns D, Ng SC, Rosenberg S, Wang X. Distinctive roles of PHAP proteins and prothymosin-alpha in a death regulatory pathway. <i>Science</i> . 2003;299(5604):223-6. Kaur R, Liu X, Gjoerup O, et al. Activation of p21-activated kinase 6 by MAP kinase kinase 6 and p38 MAP kinase. <i>J. Biol. Chem.</i> 2005; 280:3323-30.

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