

Instruction Manual

| | |
|--------------------------------|---|
| Catalog Number | PK-AB718-3149P |
| Quantity | 50 µg |
| Source | Amino acids near the amino terminus of human PHAP |
| Formulation | Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH ₂ PO ₄ , 10 mM Na ₂ HPO ₄ , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide. |
| Reconstitution | 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. |
| Storage & Stability | Store PHAP peptide at -20°C, stable for one year. |
| Application | PHAP peptide is used for blocking the activity of PHAP antibody. |
| References | 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. Alcivar A, Hu S, Tang J, et al. DEDD and DEDD2 associate with caspase-8/10 and signal cell death. <i>Oncogene</i> 2003; 22:291-7. |

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