

Instruction Manual

Catalog Number	PK-AB718-4503P
Quantity	50 µg
Source	12 amino acids near the amino terminus of human BAP3
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 BAP3 peptide at -20°C, stable for one year.
Application	BAP3 peptide is used for blocking the activity of BAP3 antibody.
References	Shiratsuchi T, Oda K, Nishimori H, et al. Cloning and characterization of BAP (BAI-associated protein 3), a C2 domain-containing protein that interacts with BAI1. <i>Biochem. Biophys. Res. Comm.</i> 1998; 251:158-65.

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