

Instruction Manual

Catalog Number	PK-AB718-2217P
Quantity	50 µg
Source	13 amino acids near the carboxy terminus of human Acinus
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 Acinus peptide at -20°C, stable for one year.
Application	Acinus peptide is used for blocking the activity of Acinus antibody.
References	Sahara S, Aoto M, Eguchi Y, Imamoto N, Yoneda Y, Tsujimoto Y. Acinus is a caspase-3-activated protein required for apoptotic chromatin condensation. <i>Nature</i> 1999 401:168-73. Geertman R, McMahon A, Sabban EL. Cloning and characterization of cDNAs for novel proteins with glutamic acid-proline dipeptide tandem repeats. <i>Biochim Biophys Acta</i> 1996;1306(2-3):147-52 (WD 0300)

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