

Instruction Manual

Catalog Number	PK-AB718-2015P
Quantity	50 µg
Source	20 amino acids near the carboxy terminus of human Apaf1
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 Apaf1 peptide at -20°C, stable for one year.
Application	Apaf1 peptide is used for blocking the activity of Apaf1 antibody.
References	Zou H, Henzel WJ, Liu X, Lutschg A, Wang X. Apaf-1, a human protein homologous to <i>C. elegans</i> CED-4, participates in cytochrome c-dependent activation of caspase-3. <i>Cell</i> 1997;90:405-13 Sakahira H, Enari M, Nagata S. Cleavage of CAD inhibitor in CAD activation and DNA degradation during apoptosis. <i>Nature</i> 1998;391:96-99 Wyllie A. Apoptosis. An endonuclease at last. <i>Nature</i> 1998;391:20-21(RD1299)

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