

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-2013P
<b>Quantity</b>	50 µg
<b>Source</b>	17 amino amino terminus of human Apaf1
<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 Apaf1 peptide at -20°C, stable for one year.
<b>Application</b>	Apaf1 peptide is used for blocking the activity of Apaf1 antibody.
<b>References</b>	<p>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.</p> <p>Alkhatib G, Liao F, Berger EA, Farber JM, Peden KW. A new SIV co-receptor, STRL33. <i>Nature</i> 1997;388:238</p> <p>Loetscher M, Amara A, Oberlin E, Brass N, Legler D, Loetscher P, D'Apuzzo M, Meese E, Rousset D, Virelizier JL, Baggiolini M, Arenzana-Seisdedos F, Moser B. TYMSTR, a putative chemokine receptor selectively expressed in activated T cells, exhibits HIV-1 coreceptor function. <i>Curr Biol</i> 1997;7:652-60 (RD1299)</p>

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