

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-1159P
<b>Quantity</b>	50 µg
<b>Source</b>	17 amino amino terminus of human FLIP
<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 FLIP peptide at -20°C, stable for one year.
<b>Application</b>	FLIP peptide is used for blocking the activity of FLIP antibody.
<b>References</b>	<p>Irmeler M, Thome M, Hahne M, Schneider P, Hofmann K, Steiner V, Bodmer JL, Schroter M, Burns K, Mattmann C, Rimoldi D, French LE, Tschopp J. Inhibition of death receptor signals by cellular FLIP. <i>Nature</i> 1997;388:190-195.</p> <p>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</p> <p>Gu J, Dong RP, Zhang C, McLaughlin DF, Wu MX, Schlossman SF. Functional interaction of DFF35 and DFF45 with caspase-activated DNA fragmentation nuclease DFF40. <i>J Biol Chem</i> 1999;274:20759-62</p>

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