

Instruction Manual

Catalog Number	PK-AB718-1161P
Quantity	50 µg
Source	18 amino acids near the carboxy terminus of human FLIP
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 FLIP peptide at -20°C, stable for one year.
Application	FLIP peptide is used for blocking the activity of FLIP antibody.
References	<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>Zhang Y, Zhou L, Miller CA. A splicing variant of a death domain protein that is regulated by a mitogen-activated kinase is a substrate for c-Jun N-terminal kinase in the human central nervous system. <i>Proc Natl Acad Sci U S A</i> 1998;95:2586-2591</p> <p>Brown TL and Howe PH. MADD is highly homologous to a Rab3 guanine-nucleotide exchange protein (Rab3-GEP). <i>Curr Biol</i> 1998;8:R191</p>

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