

Instruction Manual

Catalog Number	PK-AB718-1156
Synonyms	FLIP Antibody: I-FLICE
Description	Apoptosis is related to many diseases and induced by a family of cell death receptors and their ligands. Cell death signals are transduced by death domain (DD) containing adapter molecules and members of the ICE/CED-3 protease family. Caspases-8 (FLICE) and -10 (FLICE2) are two pivotal members in the ICE/CED-3 protease family. FLICE-inhibitory proteins were identified in virus and human and designated v-FLIPs and FLIP respectively. The human FLIP was also cloned by several labs independently and termed Casper, I-FLICE, FLAME-1, CASH, CLARP and MRIT. FLIP contains two death effector domains (DEDs) and a caspase-like domain. FLIP interacts with adapter protein FADD and caspase-8 and 10, and potently inhibits apoptosis induced by all known death receptors CD95, DR3, TRAIL-R and TNFR1.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal mFLIP antibody was raised against a peptide corresponding to amino acids 449 to 465 of mouse FLIPL/CASH (2,6).
Purification Method	Antibody is purified via DEAE-column chromatography.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Mouse
Specificity	FLIP has short form (FLIPS) and long form (FLIPL) and antibody recognizes the FLIPL only.
Formulation	Antibody is supplied in PBS containing 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	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, ICC Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. FLIP antibody can be used for the detection of FLIP by Western blot at 0.5 µg/mL. 55 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 5 µg/mL.
Images	Available upon request.
References	Hu S, Vincenz C, Ni J, Gentz R, Dixit VM. I-FLICE, a novel inhibitor of tumor necrosis factor receptor-1- and CD-95-induced apoptosis. <i>J Biol Chem</i> 1997;272:17255-17257. Thome M, Schneider P, Hofmann K, Fickenscher H, Meinl E, Neipel F, Mattmann C, Burns K, Bodmer JL, Schroter M, Scaffidi C, Krammer PH, Peter ME, Tschopp J. Viral FLICE-inhibitory proteins (FLIPs) prevent apoptosis induced by death receptors. <i>Nature</i> 1997;386:517-521. 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. Shu HB, Halpin DR, Goeddel DV. Casper is a FADD- and caspase-related inducer of apoptosis. <i>Immunity</i> 1997;6:751-763.
Images	Available upon request.
Related Products	FLIP Peptide - Cat. No. PK-AB718-1156P 3T3 (Balb) Lysate - Cat. No. PK-AB718-1212

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