

FLIP / I-FLICE (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat FLICE Inhibitory Protein

Instruction Manual

Catalog Number	PK-AB718-1161
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 and CLARP. 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. Four splice variants of c-FLIPs have been identified and termed FLIP α , β , γ , and δ , respectively.
Quantity	100 μ g
Source / Host	Rabbit
Subclass	Rabbit IgG
Immunogen	Rabbit polyclonal FLIP antibody was raised against a peptide corresponding to amino acids near the C-terminus of human FLIPaFLIPI form. The immunogen is located within the last 50 amino acids of FLIP.
Specific Species React.	Human, Mouse, Rat
Cross-Reactivity	Antibody recognizes the FLIPa only.
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Stabilizer	sodium azide
Concentration	1 mg/ml
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 μ l 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.
Purification Method	Antibody is purified via DEAE-column chromatography.
Applications	Flip antibody can be used for detection of FLIP by Western blot at 1 μ g/mL. A 55 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 10 μ g/mL. For immunofluorescence start at 10 μ g/mL.
Storage	E, WB, ICC, IF Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. 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.
References	Thome M, Schneider P, Hofmann K, Fickenscher H, Meinel 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. Nature 1997;386:517-521

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

FLIP / I-FLICE (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat FLICE Inhibitory Protein



	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. Nature 1997;388:190-195
	Shu HB, Halpin DR, Goeddel DV. Casper is a FADD- and caspase-related inducer of apoptosis. Immunity 1997;6:751-763
	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. J Biol Chem 1997;272:17255-17257
Related Products	Blocking Peptide, Cat. No. PK-AB718-1159P; K562 Lysate, Cat. No. PK-AB718-1204
	FLIP Antibody (CT), Cat. No. PK-AB718-1161; Caspase-8 Antibody (IN), Cat. No. PK-AB718-3475; Caspase-10 Antibody (CT), Cat. No. PK-AB718-1128

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