

Instruction Manual

Catalog Number	PK-AB718-2055
Synonyms	FLIP Antibody: FLIPs, FLIPg, d
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 c-FLIPs, respectively. The human FLIPs were also cloned by several labs independently and termed Casper, I-FLICE, FLAME-1, CASH, CLARP and usurpin. 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. 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 191 to 209 at C-terminus of human FLIP(/FLIPs form2.
Specific Species React.	Human
Cross-Reactivity	Antibody recognizes human and murine FLIPgamma and FLIPdelta but not FLIPalpha and FLIPbeta.
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	Affinity chromatography purified via peptide column.
Applications	FLIP can be used for detection of FLIP by Western blot at 1:1000 dilution. 35 kDa and 25 kDa bands can be detected. Antibody can also be used for immunocytochemistry starting at 2 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.

FLIPs (CT) antibody (pAb)

Rabbit Anti-Human 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	FLIP Peptide, Cat. No. PK-AB718-2055P
	Jurkat Cell Lysate, Cat. No. PK-AB718-1205

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