

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

Rabbit Anti-Human/Mouse/Rat FLICE Inhibitory Protein

Instruction Manual

Catalog Number	PK-AB718-1159
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
Immunogen	Rabbit polyclonal FLIP antibody was raised against a peptide corresponding to amino acids near the amino terminus of human FLIP. The sequence is identical in all FLIP splice variants.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	FLIP recognizes all FLIP splice variants including FLIP α , β , and γ .
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 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.
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, IHC, IF, FACS INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually. FLIP antibody can be used for Western blot at 1 - 2 μ g/mL. Immunocytochemistry at 5 μ g/mL. Antibody can also be used for immunocytochemistry starting at 5 μ g/mL. For immunofluorescence start at 20 μ g/mL.
Images	Available upon request.
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. <i>Nature</i> 1997;386:517-521 Irmler 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 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
Images	Available upon request.
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-AB718-5475; Caspase-10 Antibody (CT), Cat. No. PK-AB718-1128

FOR IN VITRO RESEARCH AND DIAGNOSTIC PROCEDURES. NOT FOR THERAPEUTIC PROCEDURES.