

Livin / KIAP antibody (pAb)

Rabbit Anti-Human Livin (Novel Member in IAP Family)

Instruction Manual

Catalog Number	PK-AB718-2505
Synonyms	Livin Antibody: KIAP
Description	Apoptosis, or programmed cell death, is related to many diseases, such as cancer. Apoptosis is triggered by a variety of stimuli including members in the TNF family and prevented by the inhibitor of apoptosis (IAP) proteins. IAP proteins form a conserved gene family that binds to and inhibits cell death proteases. A novel member in the IAP protein family was recently identified and designated Livin and KIAP for kidney IAP. Livin/KIAP contains a single baculoviral IAP repeat (BIR) domain and a RING finger domain and has two isoforms termed Livin-a and Livin-b. Transfection of Livin in cells resulted in protection from apoptosis induced by FADD, BAX, RIP, RIP3 and DR6. Livin has direct interaction with several caspases including caspase-3, -7, and -9. Livin inhibits the activation of caspase-9 induced by Apaf-1, cytochrome c, and dATP. The two isoforms of Livin appear to have different functions and tissue distributions.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Livin antibody was raised with a synthetic peptide (CRAPVRSRVRTFLS) corresponding to amino acids 264 to 280 of the short form and 281 to 298 of the long form of human Livin (1,3). This sequence is identical between (and (forms of the Livin proteins.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	A lower but much weaker band at 30 kDa was detected in Raji cell lysate, which may represent the short form of Livin.
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, IHC, IF 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.
Images	Livin antibody can be used for detection of Livin by Western blot at 0.5 µg/mL. A band at 33 kDa can be detected. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
References	Kasof GM, Gomes BC . Livin, a novel inhibitor of apoptosis protein family member. J Biol Chem. 2001;276(5):3238-46. Lin JH, Deng G, Huang Q, Morser J. KIAP, a novel member of the inhibitor of apoptosis protein family. Biochem Biophys Res Commun. 2000;279(3):820-31. Ashhab Y, Alian A, Polliack A, Panet A, Yehuda DB. Two splicing variants of a new inhibitor of apoptosis gene with different biological properties and tissue distribution pattern. FEBS Lett. 2001;495(1-2):56-60.
Images	Available upon request.
Related Products	Livin Peptide, Cat. No. PK-AB718-2505P Raji Lysat, Cat. No. PK-AB718-1207

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