

FLASH (CT) antibody (pAb)

Rabbit Anti-Human FLICE Associated Huge Protein (Homolog of CED4)

Instruction Manual

Catalog Number	PK-AB718-2271
Synonyms	FLASH Antibody: FLASH
Description	A novel mammalian CED-4 homologous was recently identified and cloned in mouse and human and designated FLASH (for FLICE-associated huge protein). FLASH is involved in Fas induced apoptosis. It is recruited to Fas after the receptor cross-linking. Overexpression of wild type of FLASH facilitates and its dominant negative form inhibits Fas induced apoptosis. FLASH interacts with the DEDs of caspase-8 and FADD through the DED-like domain of FLASH and mediates activation of caspase-8. There are parallels between FLASH and Apaf-1/CED-4 although there are arguments against their structural similarity. FLASH is widely expressed.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal FLASH antibody was raised against a synthetic peptide corresponding to amino acids near the carboxy-terminus of human FLASH (Genbank accession no. Q9UKL3), which differ from those of mouse by one amino acid.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	
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, 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. FLASH antibody can be used for detection of FLASH by Western blot at 0.5 µg/mL. An approximate 220 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 10 µg/mL. For immunofluorescence start at 10 µg/mL.
Images	Available upon request.
References	Imai Y, Kimura T, Murakami A, Yajima N, Sakamaki K, Yonehara S. The CED-4-homologous protein FLASH is involved in Fas-mediated activation of caspase-8 during apoptosis. Nature 1999;398:777-85 Koonin EV, Aravind L, Hofmann K, Tschopp J, Dixit VM Apoptosis. Searching for FLASH domains. Nature 1999;401:662-3 Medema JP. Apoptosis. Life and death in a FLASH. Nature 1999;398:756-7 (WD0800)
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2271P HeLa Cell Lysate, Cat. No. PK-AB718-1201 Apaf-1 Antibody (NT), Cat. No. PK-AB718-2013 Caspase 8 Antibody (CT), Cat. No. PK-AB718-3473

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