

# FAF1 (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat FAF1 (CT)

## Instruction Manual

Catalog Number	PK-AB718-3993
Synonyms	FAF1 Antibody: Fas-associated protein 1, hFAF1
Description	Fas-associated protein 1 (FAF1) was initially identified as a Fas-binding pro-apoptotic protein that is component of the death-inducing signaling complex in Fas-mediated apoptosis. FAF1 can also induce apoptosis in the absence of extrinsic death signals when overexpressed although it does not contain typical death motifs such as the death domain, death effector domain, and caspase recruitment domain. Overexpression of FAF1 also decreases the basal level of NF-kappaB activity in transfected 293 cells, inhibits NF-kappaB activity induced by TNF-alpha, IL-1beta and lipopolysaccharide, and prevents NF-kappaB translocation to the nucleus, suggesting that another role of FAF1 is to negatively regulate the activity of NF-kappaB. FAF1 can also interact with the inflammatory signaling PYRIN-containing Apaf-1-like proteins (PYPAFs, also called NALPs) such as PYPAF1, PYPAF2 (NALP2), and PYPAF7, suggesting FAF1 may also be involved in the inflammation pathway. Multiple differentially spliced isoforms of FAF1 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	FAF1 antibody was raised against a 19 amino acid peptide from near the carboxy terminus of human FAF1.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
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, IHC, IF Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.  FAF1 antibody can be used for detection of FAF1 by Western blot at 1 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Chu K, Niu X, and Williams LT. A Fas-associated protein factor, FAF1, potentiates Fas-mediated apoptosis. Proc. Natl. Acad. Sci. USA 1995; 92:11894-8.  Ryu SW and Kim E. Apoptosis induced by human Fas-associated factor 1, hFAF1, requires its ubiquitin homologous domain, but not the Fas-binding domain. Biochem. Biophys. Res. Commun. 2001; 286:1027-32.  Park M-Y, Jang HD, Lee SY, et al. Fas-associated Factor-1 inhibits Nuclear Factor-kappaB (NF-kappaB) activity by interfering with nuclear translocation of the RelA (p65) subunit of NF-kappaB. J. Biol. Chem. 2004; 279:2544-9.  Kinoshita T, Kondoh C, Hasegawa M, et al. Fas-associated factor 1 is a negative regulator of PYRIN-containing Apaf-1-like protein 1. Int. Immunol. 2006;18:1701-6.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-3993P; FAF1 Peptide  Cat. No. PK-AB718-1205; Jurkat Cell Lysate

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