

# Apaf-1 (NT) antibody (pAb)

## Rabbit Anti-Human/Mouse/Rat Apoptosis Protease Activating Factor 1

### Instruction Manual

Catalog Number	PK-AB718-2013
Synonyms	Apaf1 Antibody: Apaf1, Apaf-1
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 containing adapter molecules and members of the caspase family of proteases. The mammalian homologues of the key cell death gene CED-4 in <i>C. elegans</i> has been identified recently from human and mouse and designated Apaf-1 (for apoptosis protease-activating factor 1). Apaf-1 binds to cytochrome c (Apaf-2) and caspase-9 (Apaf-3), which leads to caspase-9 activation. Activated caspase-9 in turn cleaves and activates caspase-3 that is one of the key proteases, being responsible for the proteolytic cleavage of many key proteins in apoptosis. Apaf-1 can also associate with caspase-4 and caspase-8. Apaf-1 is ubiquitously expressed in human tissues.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Apaf-1 antibody was raised against a peptide corresponding to amino acids near the amino terminus of human Apaf-1. The sequences of the immunogenic peptide are identical between human and mouse.
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 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.  Apaf1 antibody can be used for detection of Apaf1 by Western blot at 1 µg/mL. A 115 - 130 kDa band should be detected. Antibody can also be used for immunohistochemistry starting at 1 µg/mL. For immunofluorescence start at 10 µg/mL.
Images	Available upon request.
References	Zou H, Henzel WJ, Liu X, Lutschg A, Wang X. Apaf-1, a human protein homologous to <i>C. elegans</i> CED-4, participates in cytochrome c-dependent activation of caspase-3. <i>Cell</i> 1997;90:405-13 Cecconi F, Alvarez-Bolado G, Meyer BI, Roth KA, Gruss P. Apaf1 (CED-4 homolog) regulates programmed cell death in mammalian development. <i>Cell</i> 1998;94:727-37 Li P, Nijhawan D, Budihardjo I, Srinivasula SM, Ahmad M, Alnemri ES, Wang X. Cytochrome c and dATP-dependent formation of Apaf-1/caspase-9 complex initiates an apoptotic protease cascade. <i>Cell</i> 1997;91:479-89 Hu Y, Benedict MA, Wu D, Inohara N, Nunez G. Bcl-XL interacts with Apaf-1 and inhibits Apaf-1-dependent caspase-9 activation. <i>Proc Natl Acad Sci USA</i> 1998;95:4386-91 (RD1299)
Images	Available upon request.
Related Products	K562 Cell Lysate, Cat. No. PK-AB718-1204 Apaf1 Antibody (CT), Cat. No. PK-AB718-2015; Caspase-9 Antibody (IN2), Cat. No. PK-AB718-2073 Caspase-4 Antibody (NT), Cat. No. PK-AB718-3451; Caspase-8 Antibody (CT), Cat. No. PK-AB718-3473

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