

PHAP (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat PHAP (Regulator of Mitochondrion Apoptosis)

Instruction Manual

Catalog Number	PK-AB718-3152
Synonyms	PHAP Antibody:
Description	Apoptosis is related to many diseases and development. Caspase-9 plays a central role in cell death induced by a variety of apoptosis activators. Cytochrome c, after released from mitochondria, binds to Apaf-1, which forms an apoptosome that in turn binds to and activate procaspase-9. Activated caspase-9 cleaves and activates the effector caspases (caspase-3, -6 and -7), which are responsible for the proteolytic cleavage of many key proteins in apoptosis. The tumor suppressor putative HLA-DR-associated proteins (PHAPs) were recently identified as important regulators of mitochondrion apoptosis. PHAP appears to facilitate apoptosome-mediated caspase-9 activation and to stimulate the mitochondrial apoptotic pathway. PHAP was also shown to oppose both Ras- and Myc-mediated cell transformation.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal PHAP antibody was raised with a synthetic peptide corresponding to amino acids at carboxy terminus of human PHAP I (GenBank Accession number P39687).
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. PHAP antibody can be used for detection of PHAP I and PHAP III by Western blot at 1 µg/mL. At approximately 35 kDa and 32 kDa a band can be detected. Antibody can also be used for immunohistochemistry starting at 10 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Jiang X, Kim HE, Shu H, Zhao Y, Zhang H, Kofron J, Donnelly J, Burns D, Ng SC, Rosenberg S, Wang X. Distinctive roles of PHAP proteins and prothymosin-α in a death regulatory pathway. Science. 2003;299(5604):223-6. Nicholson DW, Thornberry NA. Apoptosis. Life and death decisions. Science. 2003 10;299(5604):214-5.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3152P Raji Cell Lysate, Cat. No. PK-AB718-1207 Apaf-1 Antibody (CT), Cat. No. PK-AB718-2015 Caspase 9 Antibody (IN1), Cat. No. PK-AB718-2071

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