

PHAP I (IN) antibody (pAb)

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

Instruction Manual

Catalog Number	PK-AB718-3145
Synonyms	PHAP I 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 III antibody was raised with a synthetic peptide corresponding to amino acids close to carboxy terminus of human PHAP III (GenBank Accession number NP_112182).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
Specificity	This polyclonal antibody has no cross-reaction to PHAP I2a and PHAP III.
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. PHAP I antibody can be used for detection of PHAP I by Western blot at 1 - 2 µg/mL. A band at approximately 32 kDa can be detected. Antibody can also be used for immunocytochemistry starting at 2 µg/mL. For immunofluorescence start at 10 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-AB718-3145P 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.