

PHAP III antibody (pAb)

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

Instruction Manual

Catalog Number	PK-AB718-3147
Synonyms	PHAP III Antibody:
Description	Phosphoinositide 3 kinase enhancer (PIKE) is a recently identified nuclear GTPase that interacts with nuclear phosphoinositide 3-kinase (PI3 kinase) to stimulate its lipid kinase activity. PIKE exists in multiple isoforms; a shorter C-terminal isoform (PIKE-A) has also been identified as centaurin γ 1. The longest isoform (PIKE-L) has been shown to bind to the adaptor protein Homer and thereby link to metabotropic glutamate receptors, leading to activation of PI3 kinase activity and prevention of neuronal apoptosis. Overexpression of PIKE-A enhances Akt activity and promotes cancer cell invasion, whereas decreased expression of PIKE-A via dominant negative expression of PIKE-A or PIKE-A knockdown inhibits these processes. In many human cancers, expression of PIKE-A is enhanced, leading to increased Akt activity and preventing apoptosis.
Quantity	100 μ g
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, Rat
Specificity	PHAP III has no cross-reaction to PHAP I and PHAP I2a.
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 μ l 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 III antibody can be used for detection of PHAP III by Western blot at 1 μ g/mL. A band at approximately 35 kDa can be detected. Antibody can also be used for immunohistochemistry starting at 2 μ g/mL. For immunofluorescence start at 10 μ g/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-AB718-3147P A549 Cell Lysate, Cat. No. PK-AB718-1203 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.