

ZIPK (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat ZIP Kinase (Apoptosis Inducing Kinase)

Instruction Manual

Catalog Number	PK-AB718-2067
Synonyms	ZIPK Antibody: ZIPK
Description	Apoptosis is mediated by death domain containing adapter molecules and a caspase family of proteases. Certain serine/threonine protein kinases, such as ASK-1 and RIP, are mediators of apoptosis. A novel serine/threonine kinase that mediates apoptosis was recently identified and designated ZIP kinase. ZIP kinase contains an N-terminal kinase domain and a C-terminal leucine zipper structure and binds to ATF4 that is a member of ATF/CREB family. ZIP kinase has high sequence homology to DAP kinase (death-associated protein kinase), which is a mediator of apoptosis induced by γ interferon. Overexpression of ZIP kinase induces apoptosis. ZIP and DAP kinases represent a novel kinase family, which mediates apoptosis through their catalytic activities. The messenger RNA was ubiquitously expressed in various tissues.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal ZIP kinase antibody was raised against a peptide corresponding to amino acids near the center of human ZIP kinase (Genbank accession no. BAA81746).
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 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, 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. ZIPK antibody can be used for detection of DNase II expression by Western blot at 1 μ g/mL. An approximate 40 kDa band can be detected, which represents the pro-enzyme of DNase II. Antibody can also be used for immunocytochemistry starting at 10 μ g/mL. For immunofluorescence start at 10 μ g/mL.
Images	Available upon request.
References	Kawai T, Matsumoto M, Takeda K, Sanjo H, Akira S. ZIP kinase, a novel serine/threonine kinase which mediates apoptosis. Mol Cell Biol 1998;18:1642-51 (RD1299)
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2067P HeLa Lysate, Cat. No. PK-AB718-1201 ASK1 Antibody (CT), Cat. No. PK-AB718-1151 ASK1 Antibody (NT), Cat. No. PK-AB718-3679

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