

Acinus (NP) antibody (pAb)

Rabbit Anti-Human Acinus (Inducer of Chromatin Condensation)

Instruction Manual

Catalog Number	PK-AB718-2241
Synonyms	Acinus Antibody:
Description	Chromatin condensation and nuclear fragmentation (CCNF) is the hallmark of apoptosis. CCNF is triggered by the activation of members of caspase family, caspase activated DNase (CAD/DFF40), and several novel proteins including AIF and CIDE. A new inducer of chromatin condensation was recently identified and designated Acinus (for apoptotic chromatin condensation inducer in the nucleus). Acinus is cleaved by caspase-3 and an additional unknown protease generating a small active peptide p17, which causes chromatin condensation in vitro when it is added to purified nuclei. Acinus also induces apoptotic chromatin condensation in cells. Acinus is ubiquitously expressed. Three different spliced forms of Acinus have been identified in human and mouse and designated AcinusL, AcinusS and AcinusS'.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Acinus antibody was raised against a peptide (ddpvrtaqvpsprgk) corresponding to amino acids 994 to 1009 of human AcinusL, 267 to 282 of human AcinusS', or 236 to 251 of human AcinusS, which are identical to those of mouse Acinus. The selected antigenic sequence is located near the N-terminus of active peptide p17.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
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 Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. Acinus antibody can be used for detection of acinus by Western blot at 1 µg/mL. An approximate 220 kDa band can be detected.
Images	Available upon request.
References	Zamzami N, Kroemer G. Condensed matter in cell death. Nature 1999 ;401:127-8. Sahara S, Aoto M, Eguchi Y, Imamoto N, Yoneda Y, Tsujimoto Y. Acinus is a caspase-3-activated protein required for apoptotic chromatin condensation. Nature 1999 401:168-73.
Images	Available upon request.
Related Products	Acinus / NP Blocking Peptide, Cat. No. PK-AB718- 2241P K562 Lysate, Cat. No. PK-AB718-1204

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