

### Instruction Manual

|                                |  |
|--------------------------------|--|
| <b>Catalog Number</b>          | PK-AB718-3025  |
| <b>Synonyms</b>                | ARTS Antibody:   |
| <b>Description</b>             | Apoptosis is related to many diseases and development. Mitochondrial proteins, such as cytochrome c, Apaf-1, and AIF play important role in apoptosis. A novel mitochondrial septin-like protein was identified recently and designated ARTS for apoptosis related protein in TGF- $\beta$ signaling pathway. ARTS that is encoded by the human septin H5/PNUTL2/CDCrel2b gene is located to mitochondria and translocates to the nucleus when apoptosis occurs. ARTS is expressed in many tissues. It enhances cell death induced by TGF- $\beta$ and, to a lesser extent, by other apoptotic agents, such as TNF- $\alpha$ and Fas ligand.   |
| <b>Quantity</b>                | 100 $\mu$ g  |
| <b>Source / Host</b>           | Rabbit   |
| <b>Immunogen</b>               | Rabbit polyclonal ARTS antibody was raised against a synthetic peptide corresponding to amino acids near the amino terminus of human ARTS (GenBank accession number NP_536540).  |
| <b>Purification Method</b>     | Affinity chromatography purified via peptide column.   |
| <b>Clone / IgG Subtype</b>     | Polyclonal antibody  |
| <b>Species Reactivity</b>      | Human  |
| <b>Specificity</b>             |  |
| <b>Formulation</b>             | Antibody is supplied in PBS containing 0.02% sodium azide.   |
| <b>Reconstitution</b>          | During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 $\mu$ 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.   |
| <b>Storage &amp; Stability</b> | 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.   |
| <b>Applications</b>            | E, WB, IHC, IF <span style="float: right;">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.</span>   |
| <b>Images</b>                  | Available upon request.  |
| <b>References</b>              | Larisch S, Yi Y, Lotan R, Kerner H, et al. A novel mitochondrial septin-like protein, ARTS, mediates apoptosis dependent on its P-loop motif. <i>Nat Cell Biol</i> 2000;2(12):915-21<br>Depraetere V. The ARTS of apoptosis. <i>Nat Cell Biol.</i> 2000;2(12):E219.<br>Paavola,P., Horelli-Kuitunen,N., Palotie,A. and Peltonen,L. Characterization of a novel gene, PNUTL2, on human chromosome 17q22-q23 and its exclusion as the Meckel syndrome gene. <i>Genomics</i> 1999;55 (1):122-125<br>Zieger,B., Tran,H., Hainmann, I. , Wunderle,D., Zgaga-Griesz,A., Blaser,S. and Ware,J. Characterization and expression analysis of two human septin genes, PNUTL1 and PNUTL2. <i>Gene</i> 2000;261(2):197-203 |
| <b>Images</b>                  | Available upon request.  |
| <b>Related Products</b>        | Blocking Peptide, Cat. No. PK-AB718-3025P<br>Human lung Tissue Lysate, Cat. No. PK-AB718-1302<br>Apaf-1 Antibody (CT), Cat. No. PK-AB718-2015<br>AIF Antibody (CT), Cat. No. PK-AB718-2301   |

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