

Rabbit Anti-Human/Mouse ARMER (Inhibitor of Apoptosis)

Instruction Manual

Catalog Number	PK-AB718-3305
Synonyms	ARMER Antibody: ARL6IP, AIP1
Description	Apoptosis is important for normal development and tissue homeostasis. It is mediated by various caspases and ultimately results in the activation of endogenous endonucleases that degrade cellular DNA. Although apoptosis induced by endoplasmic reticulum (ER) stress is thought to be mediated by caspase-12, other caspases such as caspase-9 are also thought to be activated following ER stress. Recently, ARMER, a novel integral ER-membrane protein was shown to protect cells from ER stress-induced apoptosis. Analysis of the caspase proteolytic cascade suggests that ARMER acts by inhibiting caspase-9 activity, although the mechanism for this remains unknown. It should be noted that ARMER is not related to the inhibitor of apoptosis proteins (IAP) family and does not contain any baculoviral IAP repeat (BIR) domains.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal ARMER antibody was raised against a peptide corresponding to 15 amino acids near the C-terminus of human ARMER (GenBank accession no. Q15041).
Purification Method	Ion exchange chromatography purified
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. ARMER antibody can be used for detection of ARMER by Western blot at 0.5 to 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2 µg/mL. For immunofluorescence start at 2 µg/mL.
Images	Available upon request.
References	Stellar H. Mechanisms and genes of cellular suicide. <i>Science</i> 1995; 267:1445-9. Nakagawa T, Zhu H, Morishima N, Li E, Xu J, Yankner BA, Yuan J. Caspase-12 mediates endoplasmic-reticulum-specific apoptosis and cytotoxicity by amyloid-β. <i>Nature</i> 2000; 403:98-103. Lui HM, Chen J, Wang L, et al. ARMER, Apoptotic regulator in the membrane of the endoplasmic reticulum, a novel inhibitor of apoptosis. <i>Mol. Cancer Res.</i> 2003; 1:508-18.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3305P Mouse Small Intestine Tissue Lysate, Cat. No. PK-AB718-1408 Caspase-9 Antibody (IN1), Cat. No. PK-AB718-2071 Caspase-12 Antibody (irg), Cat. No. PK-AB718-3195
FOR IN VITRO RESEARCH ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.	