

ARC (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Apoptosis Repressor with CARD Domain

Instruction Manual

Catalog Number	PK-AB718-2185
Synonyms	ARC Antibody: ARC
Description	Apoptosis is regulated by death domain (DD) and/or caspase recruitment domain (CARD) containing molecules and a caspase family of proteases. CARD containing cell death regulators include RAIDD, RICK BCL10, Apaf-1, caspase-9, and caspase-2. A novel CARD domain containing protein was recently identified and designated ARC for apoptosis repressor with CARD . ARC interacts with caspase-2 and -8 and inhibits enzymatic activity of caspase-8. ARC suppresses apoptosis induced by cell death adapters FADD and TRADD and by cell death receptors Fas, TNFR-1, and DR3. The messenger RNA of ARC is primarily expressed in skeletal muscle and cardiac tissue .
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal ARC antibody was raised against a peptide corresponding to amino acids 2 to 18 of human origin . These sequences are identical to those of human nuclear protein Nop30 and differ from those of the rat homolog of ARC by one amino acid
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 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 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.
	ARC antibody can be used for detection of ARC by Western blot at 1:500 dilution. An approximately 25 kDa band can be detected.
Images	Available upon request.
References	Koseki T, Inohara N, Chen S, Nunez G. ARC, an inhibitor of apoptosis expressed in skeletal muscle and heart that interacts selectively with caspases. Proc Natl Acad Sci USA 1998;95:5156-60 Stoss O, Schwaiger FW, Cooper TA, Stamm S. Alternative splicing determines the intracellular localization of the novel nuclear protein Nop30 and its interaction with the splicing factor SRp30c. J Biol Chem 1999;274(16):10951-62 Geertman R, McMahon A, Sabban EL. Cloning and characterization of cDNAs for novel proteins with glutamic acid-proline dipeptide tandem repeats. Biochim Biophys Acta 1996;1306(2-3):147-52 (WD 0300)
Images	Available upon request.
Related Products	ARC Peptide, Cat. No. PK-AB718-2185P HeLa Cell Lysate, Cat. No. PK-AB718-1201

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