

PRODUCT INFORMATION



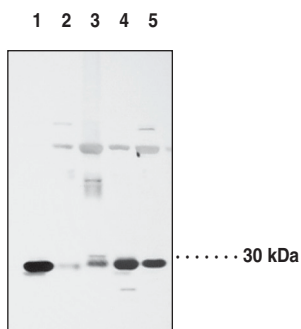
ARC Polyclonal Antibody

Item No. 160737

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
Synonym:	Apoptosis Repressor with CARD
Immunogen:	Synthetic peptide from the C-terminal region of human ARC
Species Reactivity:	(+) Human and mouse
Uniprot No.:	O60936
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 0.02% sodium azide
Host:	Rabbit
Applications:	Immunoprecipitation (IP) and Western blot (WB); the recommended starting dilution for WB is 1:200. A whole cell lysate from HeLa cells can be used as a positive control and a 25 kDa band should be detected. IP and other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Mouse cerebrum
Lane 2: Mouse liver
Lane 3: Mouse heart
Lane 4: Mouse smooth muscle
Lane 5: Mouse lung

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/20/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Apoptosis is regulated by death domain (DD) and/or caspase recruitment domain (CARD) containing molecules and a caspase family of proteases. CARD domain-containing cell death regulators included RAIDD, Apaf-1, caspase-9, and caspase-2. A novel CARD domain containing protein was recently identified and designated ARC for apoptosis repressor with CARD.¹ 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 mRNA for ARC is expressed primarily in skeletal muscle and cardiac tissues.¹

Reference

1. Koseki, T., Inohara, N., Chen, S., *et al.* ARC, an inhibitor of apoptosis expressed in skeletal muscle and heart that interacts selectively with caspases. *Proc. Nat. Acad. Sci. USA* **95**, 5156-5160 (1998).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM