

PRODUCT INFORMATION



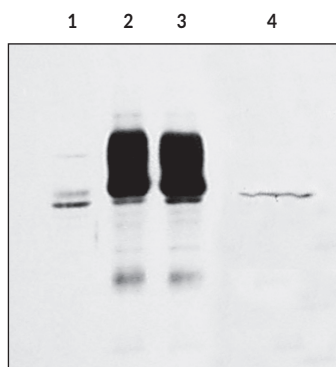
RICK Polyclonal Antibody

Item No. 160785

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
Synonyms:	CARDIAK, Rip2, Ripk2
Immunogen:	Synthetic peptide from the N-terminal region of human RICK
Species Reactivity:	(+) Human, mouse, and rat; other species not tested
Uniprot No.:	O43353
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	TBS, pH 7.4, with 50% glycerol, 0.05% BSA, and 0.02% sodium azide
Host:	Rabbit
Applications:	Western blot; the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: 293 T cell lysate
Lane 2: RICK transfected 293 T cell lysate
Lane 3: RICK transfected 293 T cell lysate
Lane 4: HeLa cell lysate

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, 11/17/2023

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 mediated by death domain (DD) and/or caspase recruitment domain (CARD) containing molecules and a caspase family of proteases. The DD-containing serine/threonine kinase RIP regulates Fas-induced apoptosis. A novel CARD-containing serine/threonine kinase that regulates apoptosis was recently identified and designated RICK for RIP-like interacting CLARP kinase.¹ RICK contains an N-terminal kinase catalytic domain and a C-terminal CARD domain. The RICK kinase domain has high sequence homology to that of RIP. Overexpression of RICK promotes the activation of caspase-8 and Fas-induced apoptosis. RICK represents a novel kinase that regulates Fas-induced apoptosis. The mRNA for RICK is expressed in multiple human tissues.¹ RICK interacts with several Nod and T-cell receptor proteins and thereby confers signal transduction for innate and adaptive immune system responses.²

References

1. Inohara, N., del Peso, L., Koseki, T., *et al.* RICK, a novel protein kinase containing a caspase recruitment domain, interacts with CLARP and regulates CD95-mediated apoptosis. *J. Biol. Chem.* **273**(29), 12296-12300 (1998).
2. Kobayashi, K., Inohara, N., Hernandez, L.D., *et al.* RICK/Rip2/CARDIAK mediates signalling for receptors of the innate and adaptive immune systems. *Nature* **416**(6877), 194-199 (2002).

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