

# RICK polyclonal antibody

The RICK (Receptor-interacting serine/threonine protein kinase 2) protein, also known as RIP-like interacting CLARP kinase, and CARD-containing ICE associated kinase, contains an N-terminal kinase domain and a C-terminal CARD domain. Overexpression of RICK results in the activation of NF- $\kappa$ B pathway and also induces apoptosis. RICK represents a kinase that regulates TNF and Fas induced-apoptosis and that is involved in the generation of proinflammatory cytokine IL-1 $\beta$ .

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 5

[View Online »](#)

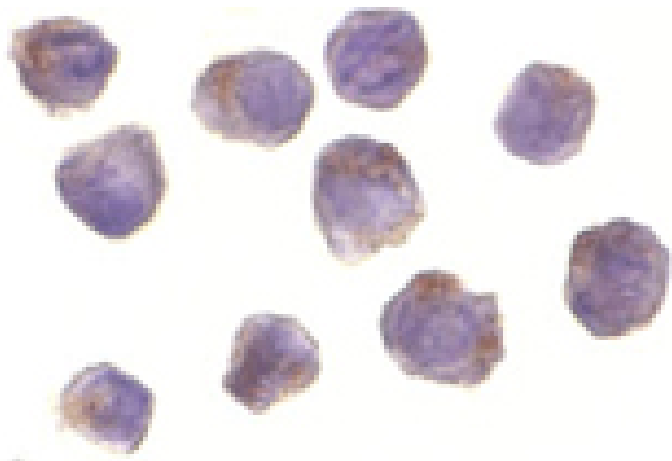
## Ordering Information

[Order Online »](#)

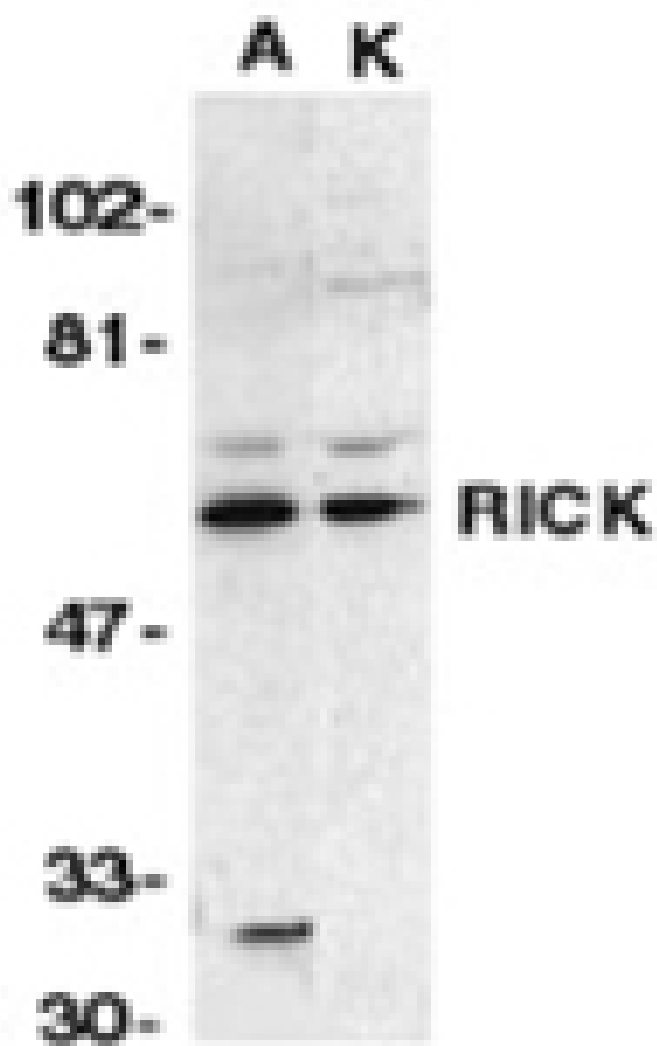
ADI-AAP-460-E	100 $\mu$ g
---------------	-------------

## Manuals, SDS & CofA

[View Online »](#)



Immunocytochemistry analysis of RICK in A431 cells with RICK antibody at 10 $\mu$ g/ml.



Western blot analysis of RICK in A431 (A) and K562 (K) whole cell lysate with RICK antibody at 1:1000 dilution.

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## Product Details

### Alternative Name

RIP2, CARDIAK, CARD-containing interleukin-1 $\beta$  converting enzyme associated kinase, RIPK2, Receptor-interacting serine/threonine protein kinase 2, Receptor-interacting protein 2

### Application

ICC, IP, WB

### Formulation

Liquid. In PBS containing 0.02% sodium azide.

### GenBank ID

AF027706

### Host

Rabbit

### Immunogen

Synthetic peptide corresponding to the sequence near the N-terminus of human RICK.

### Purity Detail

Peptide affinity purified.

### Recommendation Dilutions/Conditions

Immunocytochemistry (10 $\mu$ g/ml)Western Blot (0.5-1.0 $\mu$ g/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

### Species Reactivity

Human, Mouse, Rat

### UniProt ID

O43353

### Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)



ENZO LIFE SCIENCES,  
INC.  
May 29, 2024  
info-  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
info-  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
info-  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
info-  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
info-  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
info-  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)