

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- κ 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-1beta.

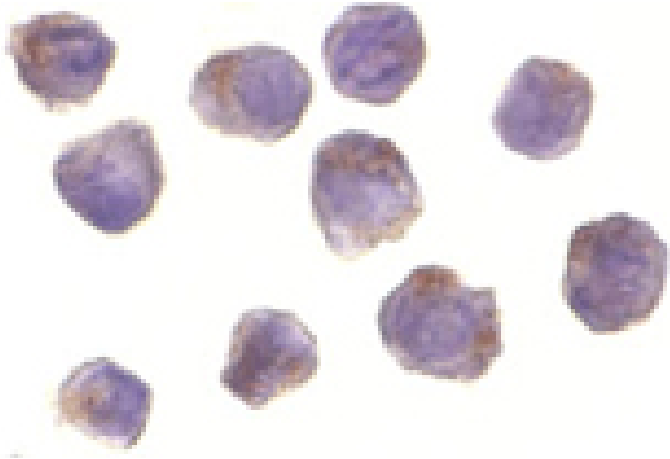
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 5

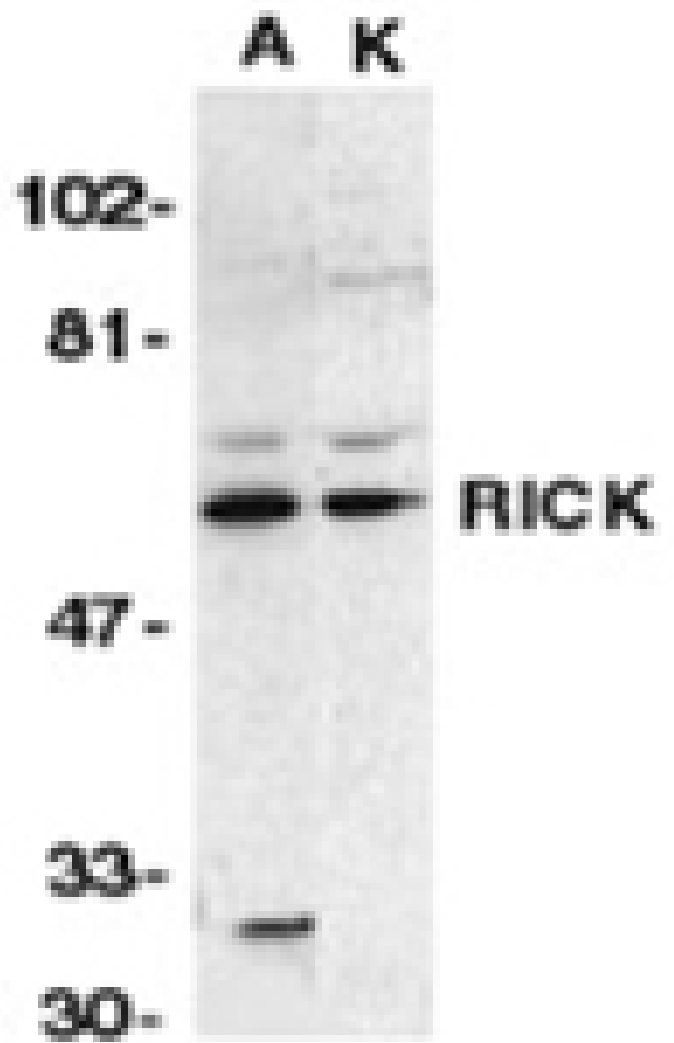
[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Immunocytochemistry analysis of RICK in A431 cells with RICK antibody at 10 μ 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 β 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 μ g/ml)Western Blot (0.5-1.0 μ 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