

# RIP3 (human) monoclonal antibody (Rippy-3)

Certain serine/threonine protein kinases, such as ASK1, RIP, DAP, and ZIP kinases, are mediators of apoptosis. Receptor interacting proteins including RIP and RIP2/RICK mediate apoptosis induced by TNF-R1 and Fas, two prototype members in the death receptor family. A member in the RIP kinase family was identified and designated RIP3. RIP3 contains an N-terminal kinase domain but, unlike RIP or RIP2, lacks the C-terminal death or CARD domain. RIP3 binds to RIP and TNF-R1, mediates TNF-R1 induced apoptosis, and attenuates RIP and TNF-R1 induced NF- $\kappa$ B activation. Overexpression of RIP3 induces apoptosis and NF- $\kappa$ B activation. The messenger RNA of RIP3 is expressed in a subset of adult tissues. Meanwhile RIP4 has been identified and initially characterized. These data suggest that RIP4 is involved in NF- $\kappa$ B and JNK signalling and that caspase-dependent processing of RIP4 may negatively regulate NF- $\kappa$ B-dependent pro-survival or pro-inflammatory signals.

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

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Receptor-interacting serine/threonine protein kinase 3, RIP-like protein kinase 3, Receptor-interacting protein 3
Application	WB
Clone	Rippy-3
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human RIP3 (receptor-interacting protein 3).
Isotype	IgG1
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human
UniProt ID	Q9Y572
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

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

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

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

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

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

Last modified: May 29, 2024