

RIP2 (human) monoclonal antibody (Nick-1)

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-139-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

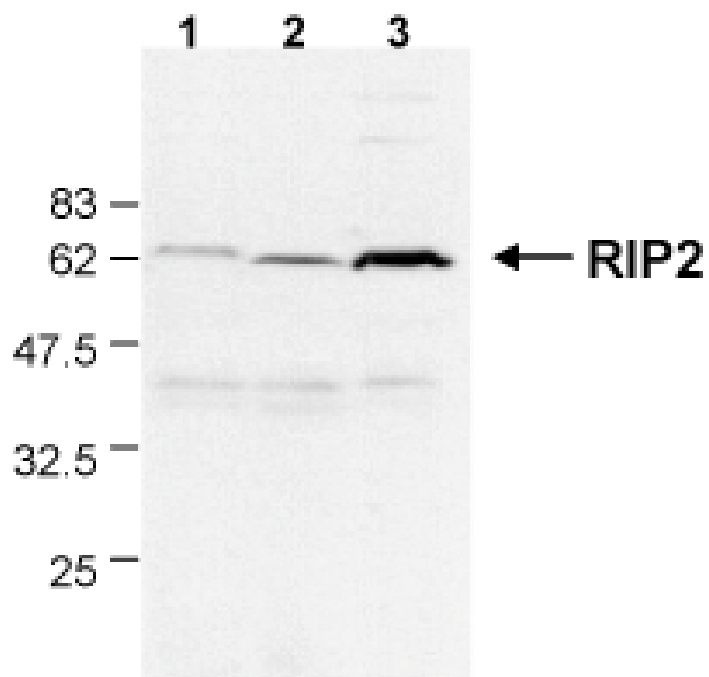


Figure: Detection of endogenous human RIP2 in different human cell lines by the MAb to RIP2 (Nick-1) (Prod. No. [ALX-804-139](#)). **Method:** Total cell extracts (50µg) from Jurkat (lane1), THP-1 (lane 2) and HT1080 (lane 3) cell lines were run on a SDS-gel, and revealed with MAb to RIP2 (Nick-1) (1:1'000 in 5% milk powder) for 1 hour at room temperature followed by incubation with HRP-coupled anti-rat secondary antibody (1:4'000).
image,RIP2 (human) monoclonal antibody (Nick-1)
image,yes

ALX-804-144,ALX-804-

144_WB.jpg,http://static.enzolifesciences.com/fileadmin/files/image/ALX-804-144_WB.jpg,Western blot analysis of ACRP30 headless (human)

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored as recommended.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	RICK, CARDIAK, CARD-containing interleukin-1 β converting enzyme associated kinase, RIPK2, Receptor-interacting serine/threonine protein kinase 2, Receptor-interacting protein 2
Application	WB
Clone	Nick-1
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant human RIP2 (receptor-interacting protein 2) (aa 1-322).
Isotype	IgG2a
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human
UniProt ID	O43353
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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

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

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

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

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

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