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 »

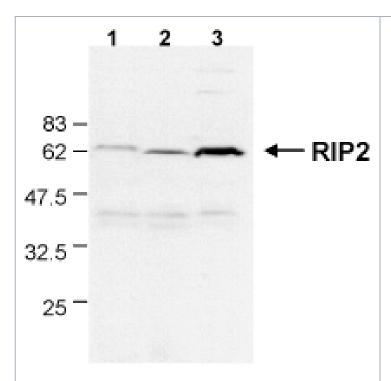


Figure: Detection of endogenous human RIP2 in different human cell lines by the MAb to RIP2 (Nick-1) (Prod. No. ALX-804-139). **Method:** Total cell extracts (50µg) from Jurkat (lane1), THP-1 (lane 2) and HT1080 (lane 3) cell lines were run on a SDS-gel, and revealed with MAb to RIP2 (Nick-1) (1:1'000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

converting enzyme associated kinase, RIPK2, Receptorinteracting 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).

lgG2a

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

UniProt ID 043353

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com