

Polyclonal Antibody to RIP3

Ordering Information

[Order Online »](#)

PSC-2283-C100	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Receptor-interacting Protein 3, Receptor-interacting Serine/Threonine Protein Kinase 3, RIP-like Protein Kinase 3

Application WB

Application Notes Detects a band of ~57kDa by Western blot.

Formulation Liquid. 100µg purified IgG in 200µl PBS containing 0.02% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 473-486 (A473QFGRGRGWQPFHK486) of mouse RIP3 (receptor-interacting protein 3).

Positive Control 3T3 whole cell lysate (Prod. No. PSC-1212) can be used as positive control.

Species Reactivity Mouse, Rat

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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