

DIY Mouse IFN Lambda 2/3 (IL-28A/B) ELISA

This DIY ELISA is a Do-It-Yourself Development kit that contains the basic components necessary for the development of sandwich ELISAs to measure natural and recombinant Human Interleukin 28A (IFN-2), and natural and recombinant human Interleukin 28B (IFN-3). VeriKine-DIY™ kits are designed for the analysis of cell culture supernates only. When used according to instructions and recommended materials, each VeriKine-DIY™ kit contains sufficient materials to develop approximately fifteen 96-well ELISA plates.

Ordering Information

[Order Online »](#)

PBL-62830-1	15x96 wells
-------------	-------------

Manuals, SDS & CofA

[View Online »](#)

- The standard in this kit is IFN Lambda 3 (IL-28B)
- This kit detects both Lambda 2 and 3 with similar reactivity

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interferon-λ 2/3, Interleukin-28A/28B

Application ELISA

Application Notes Quantitative determination of Mouse Interferon Lambda 2/3 (Interleukin-28A/B)

Assay Time Incubation time: 4 hours, 40 minutes

Contents Capture Antibody, Detection Antibody, Standard, Streptavidin-HRP

Sensitivity Assay Range: 31.25 – 2000 pg/ml

Species Reactivity Mouse

Specificity Recombinant Mouse Interferon Lambda 2 (IL-28A) and Recombinant Mouse Interferon Lambda 3 (IL-28B)



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