

DIY Human 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-61830-1	15x96 wells
-------------	-------------

Manuals, SDS & CofA

[View Online »](#)

- The standard in this kit is IFN Lambda 2 (IL-28A)
- This kit detects both Lambda 2 and 3 and is three-fold more reactive to Lambda 2 than Lambda 3

Handling & Storage

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA

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

Assay Time Incubation time: 4 hours, 45 minutes

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

Crossreactivity No cross-reactivity against Human Interferon Lambda 1 and Mouse Interferon Lambda 3.

Sensitivity Assay Range: 125 – 8000 pg/ml

Species Reactivity Human

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



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-eu@enzolifesciences.com)

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