

# DIY Human IFN Lambda 1/2/3 (IL- 29/28A/28B) 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-29 (IL-29), Human Interleukin- 28A (IL-28A), and Human Interleukin-28B (IL-28B). *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-61840-1	15x96 wells
-------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

- The standard in this kit is IFN Lambda 1 and 3
- This kit detects Lambda 1, 2 and 3, and is most reactive with Lambda 3 and 1

## 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 1/2/3 (Interleukin-29/28A/28B)

**Assay Time** Incubation time: 4 hours, 40 minutes

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

**Crossreactivity** No cross-reactivity against Human IFN- $\alpha$ A/D, IFN- $\alpha$ G, IFN- $\alpha$ A, IFN- $\alpha$ B2, IFN- $\alpha$ 1 and IFN- $\beta$ 1a.

**Sensitivity** Assay Range: 62.5 – 4000 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)