

# VeriKine-HS Human Interleukin 15 ELISA Kit

PBL's IL-15 ELISA measures both IL-15 alone and IL-15/IL15R $\alpha$  complexes in Tissue Culture Media (TCM), and complex matrices such as serum and plasma. It can also quantitate IL-15 in healthy donor matrices. The standard in this assay is the IL-15/IL-15R heterodimer but because the antibodies used are specific for IL-15, the standard curve is calibrated to the IL-15 portion of the complex. Different forms of the IL-15 receptors (IL15-R) may couple to free IL-15 protein. This assay can detect both free IL-15 and bound IL-15/IL-15R.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

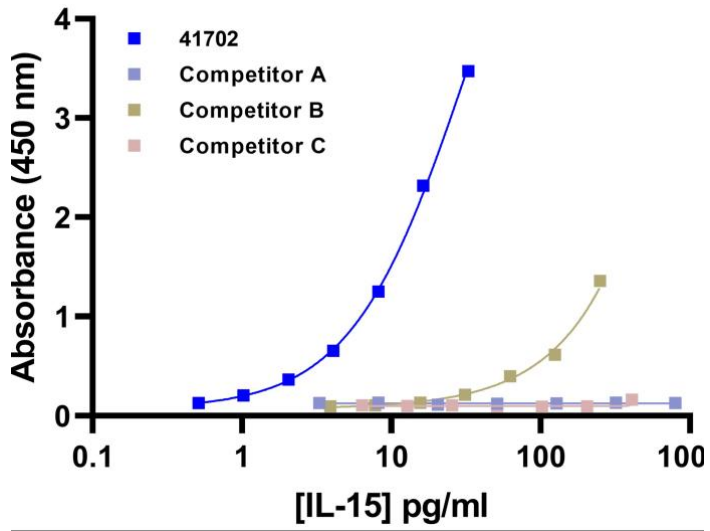
PBL-41702-1

96 wells

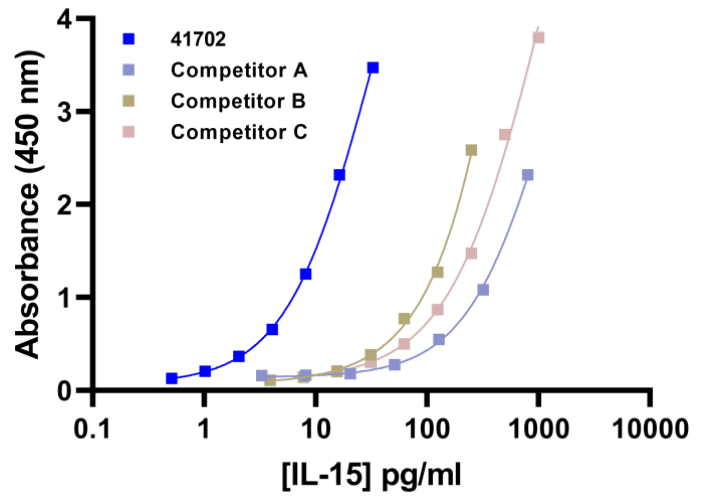
## Manuals, SDS & CofA

[View Online »](#)

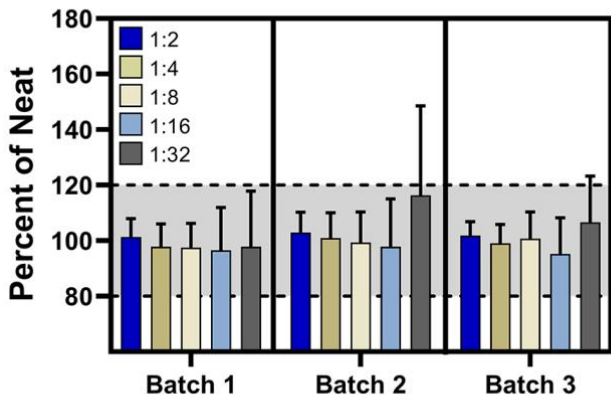
- Measures endogenous human IL-15 at or near basal level with LLOQ of 0.51 pg/ml in serum, plasma, and cell culture media
- Detects both free and bound form of IL-15 (IL-15/IL-15R complex)
- Reproducible results with inter- and intra-CVs <10% in all tested matrices



Comparison on Measuring Human IL-15/IL-15Ra complex



Comparison of Standard curves



PBL 41702 ELISA Linearity study in Normal donor serum

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name IL-15

Application ELISA

Application Notes Quantitative determination of Hu-IL-15 in serum, plasma and tissue culture media

Assay Time 4 hours

Contents Pre-Coated Micro-Titer Stripwell Plate/Sealers, Antibody Concentrate, HRP Conjugate Concentrate, Concentrate Diluent, HRP Diluent, TMB Substrate Solution, Stop Solution

Crossreactivity No cross-reactivity detected with human IL-2, IL-4, IL-7, IL-9, IL-21, IFN- $\alpha$ , IFN- $\beta$ 1a, IFN- $\gamma$ , IFN- $\omega$ ; cynomolgus IFN- $\alpha$ 2; cynomolgus/rhesus IFN- $\alpha$ ; mouse IL-15, IFN- $\alpha$ A, IFN- $\beta$ , IFN- $\gamma$ ; rat IL-15, IFN- $\alpha$ 1, IFN- $\alpha$ 14, IFN- $\beta$ , IFN- $\gamma$ ; bovine IFN-Tau2 or pig IFN- $\alpha$ .

Sensitivity 0.51 – 32.8 pg/ml

Species Reactivity Human



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)