

# VeriKine-HS Human Interleukin 22 ELISA Kit

## Ordering Information

[Order Online »](#)

PBL-41701-1	96 wells
-------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Human IL-22 ELISA with very low picogram sensitivity (LLOQ 0.78 pg/ml) for human interleukin 22 in healthy and patient donor serum, plasma, and tissue culture media (TCM)
- Accurate quantitation of endogenous human interleukin 22 in healthy and patient serum samples
- Accurate measurement of unbound (free) IL-22. IL-22 Binding Protein (IL-22BP) may interfere with the assay if present at high concentration
- Highly reproducible results in all tested matrices with inter- and intra-CVs < 10%

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name IL-22

Application ELISA

Application Notes Quantitative determination of Hu-IL-22 in serum, plasma and tissue culture media

Assay Time 4 hours

Contents Pre-Coated Micro-Titer Stripwell Plate/Sealers, Wash Solution Concentrate, Human Interleukin-22 Standard (10,000 pg/ml), Standard Diluent, Assay Buffer, Antibody Concentrate, HRP Conjugate Concentrate, Concentrate Diluent, HRP Diluent, TMB Substrate Solution, Stop Solution

Crossreactivity Cross-reacts with Monkey IL-22 (< 0.03%). No cross-reactivity detected with human IL-10, IL-17, IL-19, IL-20, IL-24, IL-26/AK155, IFN-β1a, IFN-γ, IFN-ω; cynomolgus IFN-α2; cynomolgus/rhesus IFN-α; mouse IL-22, IFN-αA, IFN-β, IFN-γ; rat IL-22, IFN-α1, IFN-α14, IFN-β, IFN-γ; bovine IFN-Tau2 or pig IFN-α.

Sensitivity 0.78 – 50 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-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-de@enzolifesciences.com](mailto:info-de@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)