

# VeriKine Human Interferon Alpha ELISA Kit

## Quantifies human IFN-alpha in cell or tissue culture media down to 12.5 pg/ml

The VeriKine™ Human Interferon Alpha ELISA Kit has been developed to quantify levels of human IFN-alpha in cell or tissue culture media via a sandwich enzyme-linked immunoassay format.

Citations: 3

[View Online »](#)

## Ordering Information

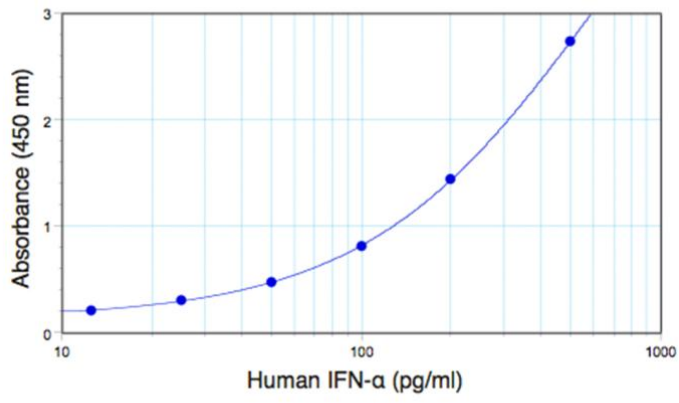
[Order Online »](#)

PBL-41100-1	96 wells
PBL-41100-2	5x96 wells

## Manuals, SDS & CofA

[View Online »](#)

- ELISA assay quantifies human IFN alpha in tissue culture media (TCM) down to 12.5 pg/ml
- Less than 10% inter- and intra-assay CV
- Less than 4 hours assay time



Representative IFN-Alpha standard curve from 12.5 pg/ml to 500 pg/ml

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name IFN- $\alpha$ , Interferon- $\alpha$

Application ELISA

Application Notes Quantitative determination of Hu-IFN alpha A, 2, A/D, D, K and 4b in tissue culture media

Assay Time Incubation time: 3 hours 15 minutes

Contents Pre-coated microtiter plate(s), Plate Sealers, Wash Solution Concentrate, Human Interferon Alpha Standard (10,000 pg/ml), Dilution Buffer, Antibody Concentrate, HRP Conjugate Concentrate, Concentrate Diluent, TMB Substrate, Stop Solution

Crossreactivity No cross-reactivity against human IFN-beta, gamma, omega; rat IFN-alpha, beta, gamma; mouse IFN-alpha, beta, gamma; or bovine IFN-tau.

Sensitivity 12.5 – 1000 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)