

# VeriKine Human Interferon Alpha Multi-Subtype Serum ELISA Kit

The VeriKine™ Human Interferon Alpha Multi-Subtype Serum ELISA kit is designed to provide accurate detection of human IFN-alpha in serum and plasma samples.

Citations: 4

[View Online »](#)

## Ordering Information

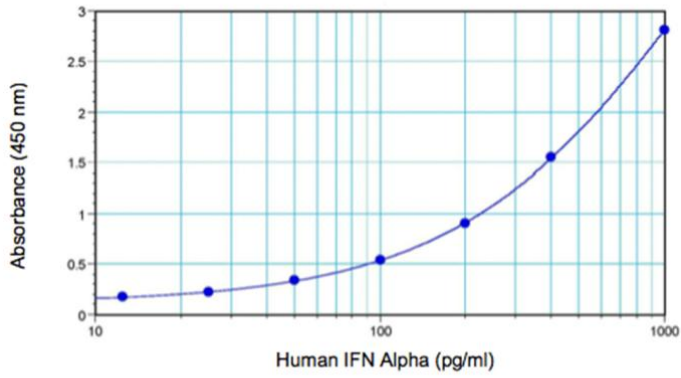
[Order Online »](#)

PBL-41110-1	96 wells
PBL-41110-2	96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Human IFN-Alpha ELISA kit for measuring global human interferon alpha amount in serum, plasma, and tissue culture media (TCM)
- Greater than 90% spike recovery in normal human serum
- ELISA assay detects 11 of 12 major human IFN-alpha subtypes
- Less than 10% inter and intra-assay CV



Representative IFN-Alpha standard curve from 12.5 pg/ml to 1,000 pg/ml in standard diluent

Spike Recovery Matrix	Low (30 pg/ml)	Medium (300 pg/ml)	High (800 pg/ml)
Normal human serum	94%	100%	96%
Normal human plasma (Na-Citrate)	70%	112%	110%
Normal human plasma (Na-Heparin)	95%	106%	101%
Normal human plasma (Na-EDTA)	60%	86%	84%

Spike-Recovery of IFN-Alpha in different matrices

## 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, B2, C, D, G, H, I, J1, K, 1, 4b and WA 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), Sample Diluent, Antibody Concentrate, HRP Conjugate Concentrate, Concentrate Diluent, TMB Substrate, Stop Solution, Dilution Buffer

Crossreactivity No cross reactivity with human interferon beta, gamma, omega; mouse or rat interferon alpha, beta, gamma; or bovine interferon 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)