

VeriKine-HS Human Interferon Alpha All Subtype ELISA Kit

This high sensitivity Human IFN Alpha ELISA kit is designed to provide accurate low pg/ml measurement of ALL human interferon alpha subtypes in autoimmune sera, normal serum, EDTA plasma, and tissue culture media (TCM).

Citations: 5

[View Online »](#)

Ordering Information

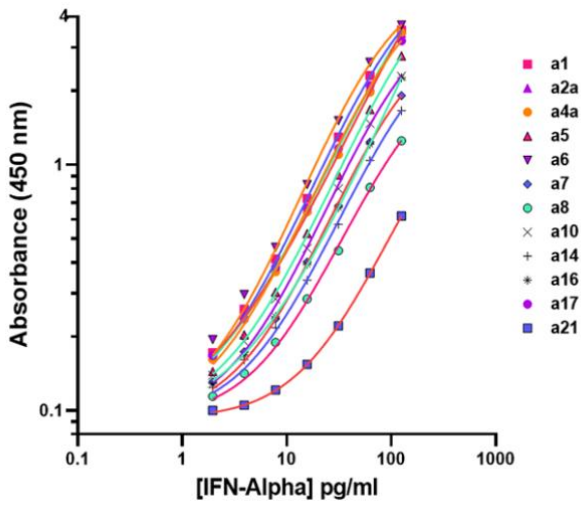
[Order Online »](#)

PBL-41115-1	96 wells
-------------	----------

Manuals, SDS & CofA

[View Online »](#)

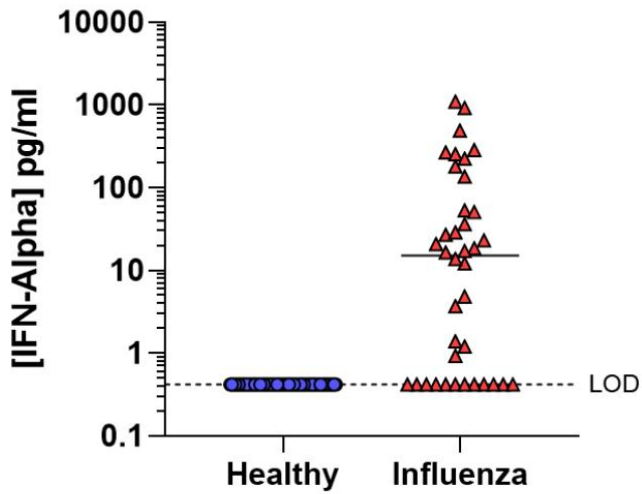
- Human IFN Alpha ELISA quantitates low level human interferon alpha in normal & autoimmune sera, EDTA plasma, and TCM with LLOQ 1.95 pg/ml
- Detects All 12 human IFN Alpha subtypes for accurate total IFN alpha measurement in samples
- Reproducible inter- (<10%) and intra-assay (< 8%) CVs in all tested matrices provide reliable data



Representative Human IFN-Alpha Subtype standard curves in Healthy Donor Serum (HDS)

Subtype	LOD (pg/ml)	LOQ (pg/ml)
$\alpha 1$ (αD)	0.09	0.343
$\alpha 2a$ (αA)	0.056	0.214
$\alpha 4a$ ($\alpha M1$)	0.099	0.382
$\alpha 5$ (αG)	0.083	0.276
$\alpha 6$ (αK)	0.041	0.183
$\alpha 7$ ($\alpha J1$)	0.112	0.423
$\alpha 8$ ($\alpha B2$)	0.284	0.993
$\alpha 10$ (αC)	0.108	0.447
$\alpha 14$ (αH)	0.168	0.585
$\alpha 16$ (αWA)	0.078	0.328
$\alpha 17$ (αI)	0.156	0.522
$\alpha 21$ (αF)	0.293	1.129

Representative Lower Limits of Quantification (LLOQ) and Lower Limits of Detection (LLOD) in Healthy Donor Serum for each human IFN-Alpha Subtype



IFN-Alpha level in Healthy and Influenza patient serum samples measured by PBL 41115-1 ELISA

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IFN- α , Interferon- α

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 serum, plasma and tissue culture media

Assay Time Incubation time: 22 hours, 30 minutes

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

Crossreactivity Cross-reacts with Cynomolgus/Rhesus IFN- α . No cross-reactivity against human IFN- β , IFN- γ , or IFN- ω ; mouse or rat IFN- α , IFN- β or IFN- γ ; or bovine IFN- τ .

Sensitivity 1.95-125 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)