

VeriKine Human Interferon Alpha Multi-Subtype ELISA Kit

Detects 14 out of 15 human interferon alpha subtypes with global IFN alpha measurement

The VeriKine Human IFN- α Multi-Subtype ELISA kit is specifically formulated to detect 14 out of 15 identified human IFN- α subtypes/alleles in media using a sandwich immunoassay.

Ordering Information

[Order Online »](#)

PBL-41105-1	96 wells
PBL-41105-2	5x96 wells

Manuals, SDS & CofA

[View Online »](#)

- Human IFN Alpha ELISA assay detects 11 of 12 human IFN alpha subtypes
- Assay Ranges from 12.5 - 500 pg/ml and from 156 - 5000 pg/ml
- Less than 10% inter and intra-assay CV

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 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 -500 pg/ml / 156 – 5000 pg/ml

Species Reactivity Human



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com