

VeriKine-HS Human Interferon Beta Serum ELISA Kit

Detects Human IFN beta in serum, plasma or cell supernatant as low as 2.3 pg/ml

The VeriKine-HS™ Human Interferon Beta Serum ELISA Kit has been developed to measure low/basal levels of human IFN-β in tissue culture media, plasma or serum samples by sandwich enzyme linked immunosorbent assay (ELISA).

Citations: 5

[View Online »](#)

Ordering Information

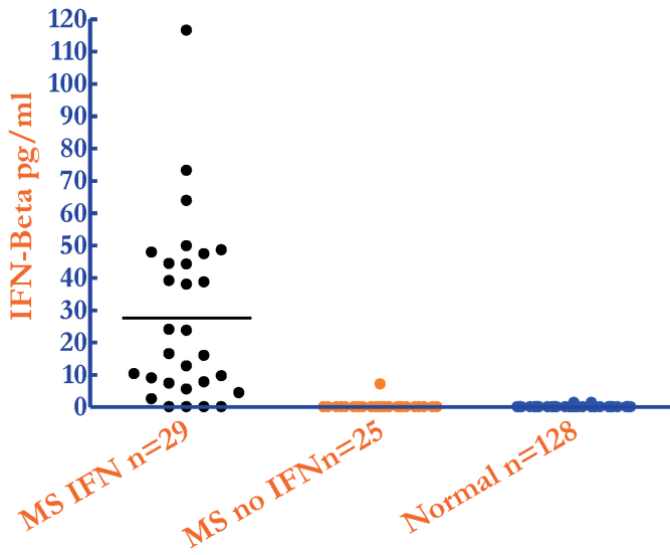
[Order Online »](#)

PBL-41415-1	96 wells
-------------	----------

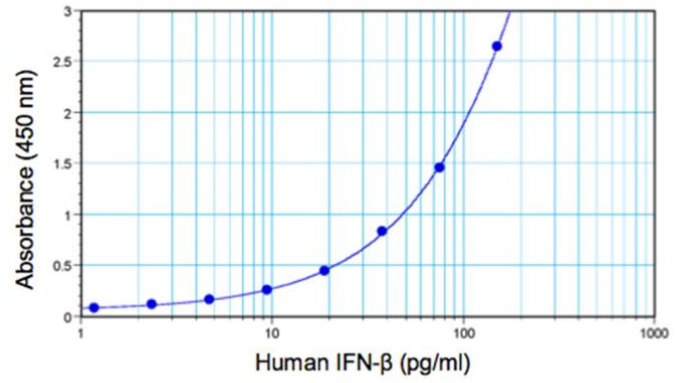
Manuals, SDS & CofA

[View Online »](#)

- IFN-beta ELISA is designed to measure human interferon beta in serum, plasma, and tissue culture media with 1.2 pg/ml LLOQ sensitivity
- Suitable for measurement of IFN-beta in autoimmune serum
- Suitable for measurement of trademarked interferon beta 1a and interferon beta 1b therapeutic molecules in human serum
- ELISA assay provides > 90% serum spike-recovery
- Reproducible results with inter-assay CV < 8% and intra-assay CV < 10%
- No inhibition of IFN beta detection by 3-log excess of soluble IFNAR2 protein



Levels of IFN-Beta in 128 Normal donor, 54 Multiple Sclerosis (MS) patient sera on and off IFN-Beta therapies. IFN-Beta is quantified in 865 of MS patients on IFN-Beta therapy, 4.1% of MS patients on other therapies and 1.5% of Normal donor sample



Representative Human IFN-Beta standard curve from 1.2 – 150 pg/ml

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IFN- β , Interferon- β

Application ELISA

Application Notes Quantitative determination of Hu-IFN Beta in tissue culture media, serum and plasma

Assay Time Incubation time: 3 hours 30 minutes

Contents Pre-coated microtiter plate, Plate Sealers, Wash Solution Concentrate, Human Interferon Beta 1a Standard (100,000 pg/ml), Standard Diluent, Sample Buffer, Antibody Concentrate, HRP Conjugate Concentrate, Diluent Additive III, Assay Diluent, TMB Substrate, Stop Solution

Crossreactivity No cross reactivity against human IFN Alpha, IFN Gamma, IFN Omega, IL-6, mouse IFN Alpha, IFN Beta or rat IFN Beta.

Sensitivity 1.2 – 150 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