

VeriKine Human Interferon Beta ELISA Kit

Quantifies human IFN Beta in tissue culture media to as low as 25 pg/ml

The VeriKine™ Human Interferon Beta ELISA Kit has been developed to measure levels of human IFN- β in tissue culture media samples by sandwich enzyme linked immunosorbent assay (ELISA).

Citations: 5

[View Online »](#)

Ordering Information

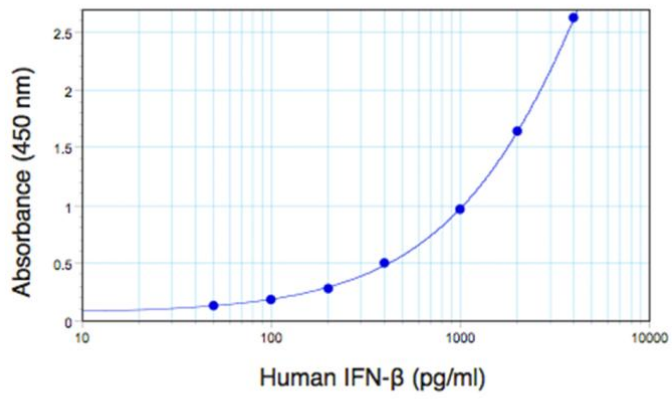
[Order Online »](#)

PBL-41410-1	96 wells
PBL-41410-2	5x96 wells

Manuals, SDS & CofA

[View Online »](#)

- Quantifies human IFN beta in tissue culture media (TCM)
- Reproducible results with inter-assay CV < 8% and intra-assay CV < 8%
- Spike recovery > 90%
- Recombinant Human IFN beta 1a expressed in mammalian cells provided as the standard



Typical Human IFN-Beta standard curve from 50 to 4,000 pg/ml using PBL's Human IFN-Beta ELISA (41410)

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

Assay Time Incubation time: 3 hours 15 minutes

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

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

Sensitivity 50 – 4000 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