

VeriKine-HS Human Interferon Beta TCM ELISA Kit

The VeriKine-HS Human IFN-Beta TCM ELISA Kit is designed for accurate quantification of low level human interferon beta in buffers or tissue culture media (TCM) by sandwich ELISA.

Citations: 4

[View Online »](#)

Ordering Information

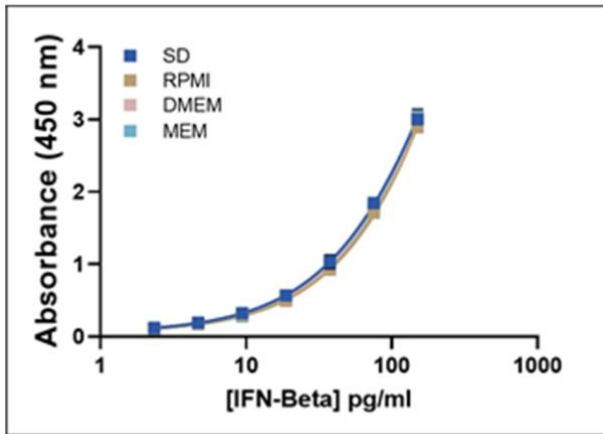
[Order Online »](#)

PBL-41435-1	96 wells
-------------	----------

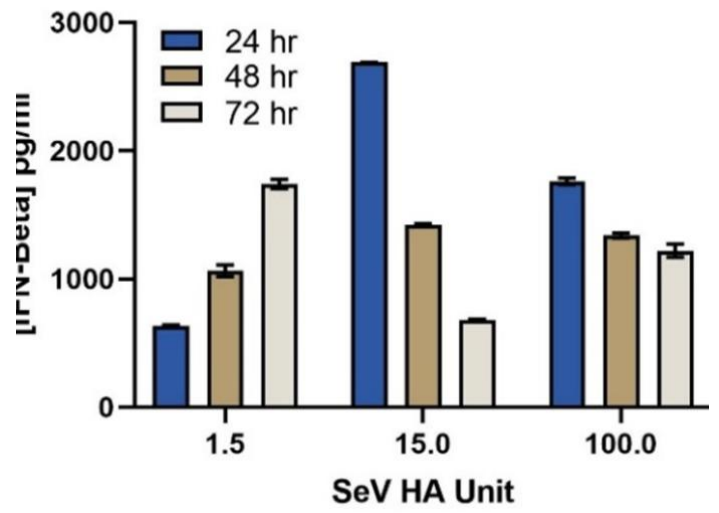
Manuals, SDS & CofA

[View Online »](#)

- Human IFN-Beta ELISA detects the amount of interferon beta in tissue culture media (TCM) to as low as 2.34 pg/ml
- Total IFN-Beta quantification with no interference from soluble IFN receptor (sIFNAR2)
- Reproducible results with inter- and intra-assay CVs < 10%



Representative Human IFN-Beta standard curve in various media matrices



Endogenous levels of Human IFN-Beta quantified in Sendai virus induced HEK293 Cells

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 High sensitivity detection of Hu IFN-Beta in tissue culture media

Assay Time 3 hours

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

Crossreactivity No cross-reactivity with human IFN-alpha, IFN-gamma, IFN-lambda, IL-6; mouse IFN-alpha, IFN-beta; or rat IFN-beta

Sensitivity 2.34 – 150 pg/ml

Species Reactivity Human

Specificity Human IFN Beta 1a & 1b.



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