

VeriKine-HS Mouse Interferon Beta Serum ELISA Kit

The world's most sensitive Mouse IFN Beta ELISA

PBL's 96-well high sensitivity ELISA kit will equip you with the most sensitive and consistent tool for accurately measuring Mouse IFN- β in serum and plasma matrices. This kit quantitates Mouse Interferon Beta (IFN- β) in sera, plasma and tissue culture media by sandwich enzyme linked immunosorbent assay (ELISA). Interferon binds to plates coated with antibody and detection is accomplished using a detection antibody followed by streptavidin conjugated to horseradish peroxidase (HRP). This ELISA kit utilizes Tetramethyl-benzidine (TMB) as the substrate. Purified, recombinant mouse IFN- β expressed in mammalian cells is provided as the standard. The sensitivity provided by this kit, to less than 1 pg/ml in serum and plasma, will propel your mouse studies and help uncover insights into the pathology of diseases.

Citations: 5

[View Online »](#)

Ordering Information

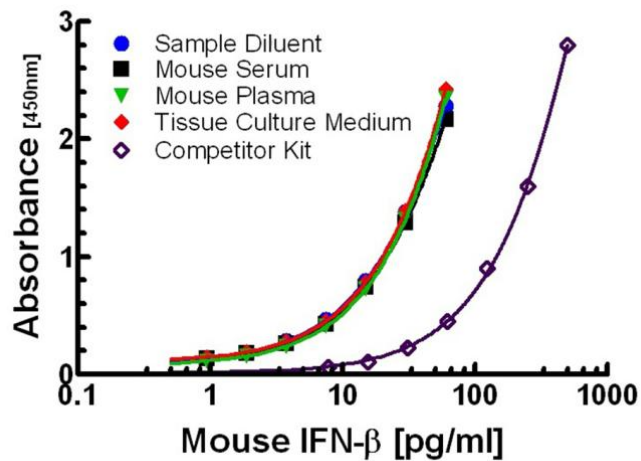
[Order Online »](#)

PBL-42410-1	96 wells
PBL-42410-2	5x96 wells

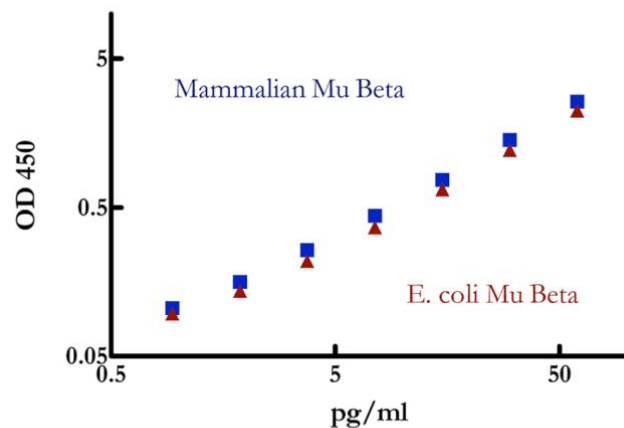
Manuals, SDS & CofA

[View Online »](#)

- Mouse IFN-Beta ELISA with 0.94 pg/ml LLOQ for accurate quantitation of mouse interferon beta protein
- Compatible with serum, plasma, and tissue culture media (TCM)
- Provides reproducible results with inter- & intra-assay CVs of < 10%
- Total assay time of ~2 hours



Typical standard curve in a variety of matrices for PBL's High-Sensitivity Mouse IFN-Beta ELISA (42410) and another supplier's ELISA in assay buffer



Comparison of mammalian Cell- and E.coli-expressed Mouse IFN-Beta detection using PBL's 42410 ELISA kit

Concentration (pg/ml)	2.5	10	50
Intra-Assay CV	5.7%	5.5%	5.5%
Inter-Assay CV	8.3%	6.3%	4.8%
Average Recovery	98%	92%	94%

Intra- and Inter-assay CVs and average percent recoveries using PBL's high sensitivity Mouse IFN-Beta ELISA (42410)

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 Mu-IFN Beta in serum, plasma and tissue culture media

Assay Time Incubation time: 1 hour, 50 minutes

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

Crossreactivity No cross-reactivity against mouse: IFN- α A, IFN- α 4, IFN- α 11, IFN- α 13, IFN- γ , IL-6, Limitin, TNF- α ; human: IFN- α A, IFN- β , IFN- γ , IFN- ω ; rat: IFN- α , IFN- β , IFN- γ ; swine: IFN- α , IFN- β

Sensitivity 0.94 – 60 pg/ml

Species Reactivity Mouse



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