

# VeriKine Mouse Interferon Beta ELISA Kit

## Quantifies mouse interferon beta in tissue culture or serum to as low as 15.6 pg/ml

The VeriKine™ Mouse Interferon Beta ELISA Kit has been developed to quantitate levels of IFN- $\beta$  in tissue culture media or serum samples in a sandwich enzyme immunoassay format with recombinant mouse IFN-beta expressed in mammalian cells provided as the standard in the kit.

Citations: 5

[View Online »](#)

## Ordering Information

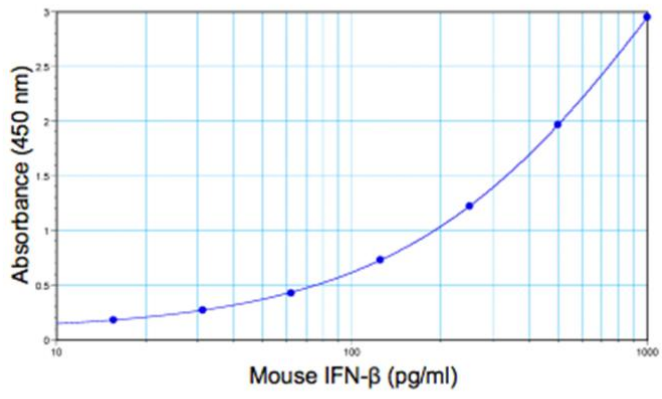
[Order Online »](#)

|             |            |
|-------------|------------|
| PBL-42400-1 | 96 wells   |
| PBL-42400-2 | 5x96 wells |

## Manuals, SDS & CofA

[View Online »](#)

- Measures mouse IFN beta in tissue culture media (TCM) and mouse serum
- Less than 10% intra- and Inter-assay CVs



Representative standard curve from 15.6 to 1,000 pg/ml for PBL's Mouse IFN-Beta ELISA (42400)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name IFN- $\beta$ , Interferon- $\beta$

Application ELISA

Application Notes Quantitative determination of Mu-IFN Beta in tissue culture media and serum

Assay Time Incubation time: 3 hours 15 minutes

Contents Pre-coated microtiter plate, Plate Sealers, Wash Solution Concentrate, Mouse Interferon Beta Standard (500,000 pg/ml), Dilution Buffer, Antibody Concentrate, HRP Conjugate Concentrate, Concentrate Diluent, TMB Substrate, Stop Solution

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

Sensitivity 15.6 – 1000 pg/ml

Species Reactivity Mouse



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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