

# VeriKine Mouse Interferon Alpha ELISA Kit

**Detects global amount of mouse interferon alpha in tissue culture media as low as 12.5 pg/ml**

The VeriKine™ Mouse Interferon Alpha ELISA Kit has been developed to quantitate levels of IFN-alpha in tissue culture media using a sandwich immunoassay.

Citations: 5

[View Online »](#)

**Ordering Information**

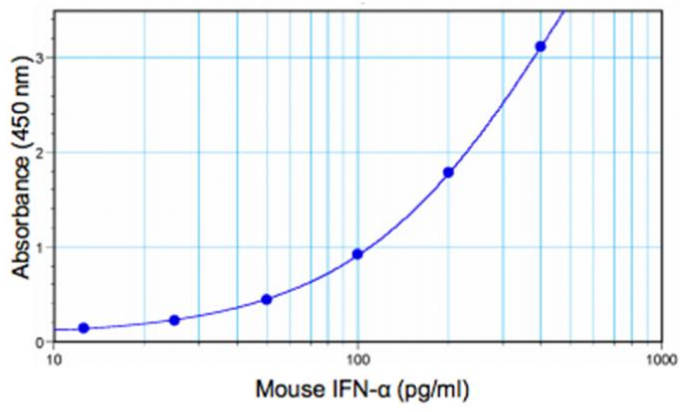
[Order Online »](#)

PBL-42120-1	96 wells
PBL-42120-2	5x96 wells

**Manuals, SDS & CofA**

[View Online »](#)

- Detects all 14 mouse IFN Alpha subtypes in tissue culture media (TCM)
- Consistent performance and reproducible results with < 10% CVs



Typical Mouse IFN Alpha standard curve from 12.5/ml to 400 pg/ml

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

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

Application ELISA

Application Notes Quantitative determination of Mu-IFN Alpha in tissue culture media

Assay Time Incubation time: 27 hours, 15 minutes

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

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

Sensitivity 12.5 – 400 pg/ml

Species Reactivity Mouse

Specificity Detects all 14 mouse IFN-alpha sub-types in tissue culture media.



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)