

VeriKine-HS Mouse Interferon Alpha All Subtype ELISA Kit

The VeriKine-HS Mouse IFN- α ELISA kit has been developed to quantitate low levels of IFN- α in mouse serum, plasma and cell/tissue culture media (10% FBS) using a sandwich immunoassay. The kit contains an anti-mouse IFN- α detection antibody conjugated to biotin and quantified by streptavidin horseradish peroxidase (HRP) using tetramethyl-benzidine (TMB) as a substrate. The kit standard is mouse IFN- α 4 protein, and typical standard curves are included with the procedure. It is important to run the standard curve in the same matrix as one's samples unless one has demonstrated that the matrix does not significantly affect the signal.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

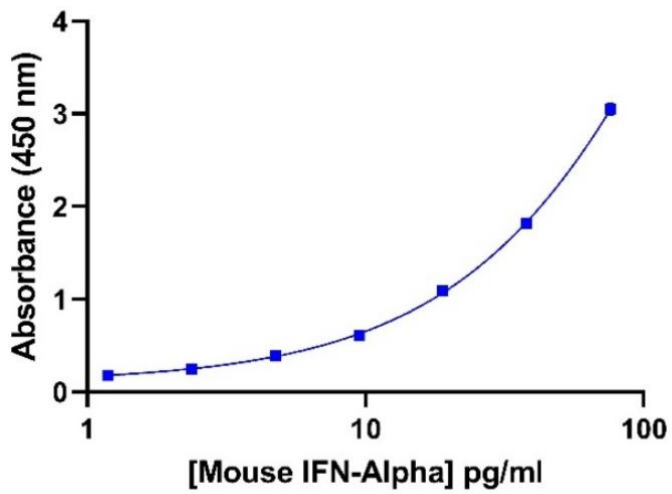
PBL-42115-1

96 wells

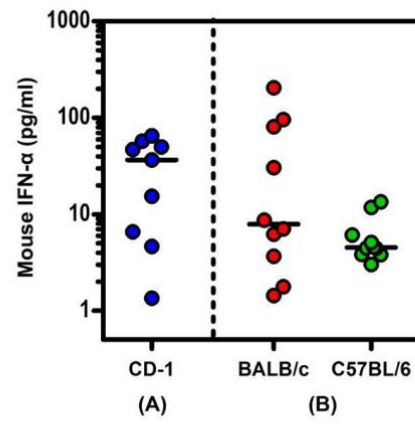
Manuals, SDS & CofA

[View Online »](#)

- Sensitive global detection of **all 14** mouse IFN alpha in **serum, plasma**, and tissue culture media
- Quantification of mouse IFN- α as low as **2.38 pg/ml**
- Rapid ELISA provides total interferon alpha readout in **~ 2 hrs**
- High data reproducibility with **less than 10% CVs**



Representative Mouse IFN-Alpha standard curve in sample diluent



Levels of endogenous mouse IFN-α quantified in pooled plasma (A) and individual sera (B) using 42115-1 ELISA

Endogenous Mouse IFN-Alpha Quantified in plasma and serum from three popular mouse strains using PBL's high sensitivity Mouse IFN-Alpha All-Subtype ELISA (42115)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IFN-alpha

Application ELISA

Application Notes Quantitative determination of Mu-IFN Alpha in serum, plasma, and tissue culture media

Assay Time Typical Incubation time: 1 hour, 54 minutes

Contents Pre-coated microtiter plate, Plate Sealers, Wash Solution Concentrate, Mouse Interferon Alpha 4 Standard (10,000 pg/ml), Sample Diluent, Antibody Concentrate, Antibody Diluent, HRP Conjugate Concentrate, HRP Diluent, TMB Substrate Solution, Stop Solution

Crossreactivity No cross-reactivity against mouse IFN- β , mouse IFN- γ , rat IFN- α 1, or IFN- α 14; very weak cross-reactivity with human IFN- α 14.

Sensitivity 2.38 – 152 pg/ml

Species Reactivity Mouse

Specificity Recognizes all 14 Mouse Interferon Alpha subtypes.



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