

VeriKine

Cynomolgus/Rhesus

Interferon Alpha

Serum ELISA Kit

The Cynomolgus/Rhesus ELISA Kit has been developed to analyze the presence of non-human primate IFN-alpha in tissue culture media or serum samples by sandwich enzyme linked immunosorbent assay (ELISA).
Quantifies non-human primate IFN alpha in serum, plasma, or tissue culture media to as low as 25 pg/ml

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

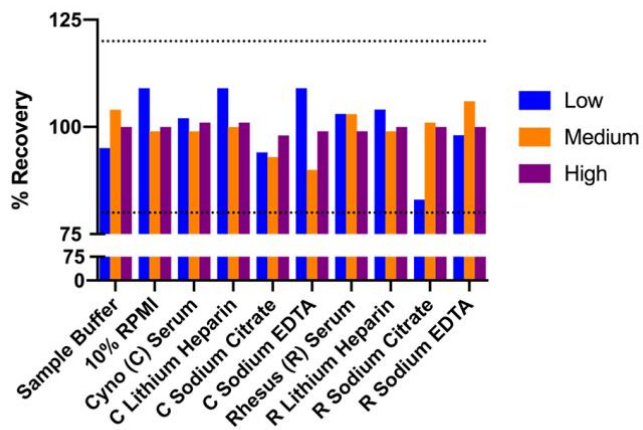
PBL-46100-1

96 wells

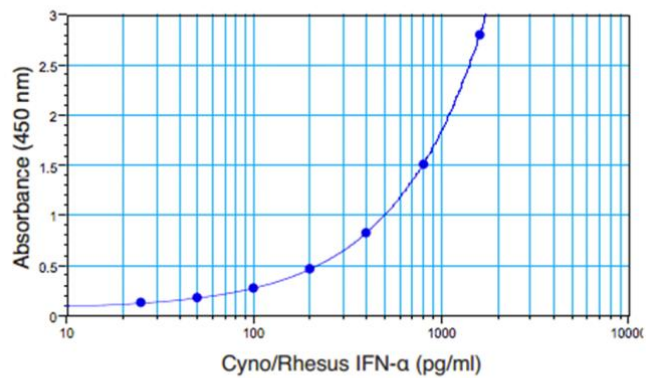
Manuals, SDS & CofA

[View Online »](#)

- Quantifies Cynomolgus and Rhesus IFN-Alpha 2 in serum, plasma, and tissue culture media (TCM)
- Reproducible results with tissue culture media inter-assay CV \leq 9.2% and intra-assay CV \leq 2.5% (serum \leq 14.4% and \leq 9.2%, respectively)
- Recombinant Rhesus/Cyno IFN Alpha 2 provided as the ELISA standard



Spike-recovery of Cynomolgus/Rhesus IFN-Alpha in serum, plasma, TCM and 46100 kit's sample buffer



Representative Cynomolgus IFN-Alpha standard curve from 25-1,600 pg/ml

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IFN- α , Interferon- α

Application ELISA

Application Notes Quantitative determination of Cynomolgus/Rhesus Macaque IFN Alpha in tissue culture media, serum and plasma

Assay Time Incubation time: 3 hours 15 minutes

Contents Pre-coated microtiter plate, Plate sealers, Wash Solution Concentrate, Rhesus/Cyno Interferon Alpha 2 Standard (10,000 pg/ml), Standard Diluent, Sample Buffer, Antibody Concentrate, HRP Conjugate Concentrate, Concentrate Diluent, TMB Substrate, Stop Solution

Crossreactivity Strong cross-reactivity against human IFN alpha2; No cross-reactivity with human IFN gamma; or mouse and rat IFN alpha.

Sensitivity 25 – 1600 pg/ml

Species Reactivity Cynomolgus, Rhesus monkey



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