

VeriKine Cynomolgus Interferon Beta ELISA Kit

This kit has been developed to measure low/basal levels of *Macaca fascicularis* IFN- β in a variety of sample matrices including serum, plasma and tissue culture media.

Ordering Information

[Order Online »](#)

PBL-46415-1	96 wells
-------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Quantifies Cynomolgus IFN-Beta in serum, plasma, and tissue culture media (TCM) as low as 5.47 pg/ml
- Reproducible results with inter- and intra-assay CVs \leq 10%
- Recombinant Mammalian Cynomolgus IFN Beta provided as the ELISA standard

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 Cynomolgus IFN Beta in tissue culture media, serum and plasma

Assay Time Incubation time: 3 hours

Contents Pre-coated microtiter plate, Plate Sealers, Wash Solution Concentrate, Cynomolgus Interferon Beta Standard (100,000 pg/ml), Standard Diluent, Sample Buffer, Antibody Concentrate, HRP Conjugate Concentrate, Assay Diluent, TMB Substrate Solution, Stop Solution

Sensitivity 5.47 – 350 pg/ml

Species Reactivity Cynomolgus



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