

# VeriKine Pig Interferon Alpha Subtype ELISA Kit

The Verikine-HS Pig IFN- $\alpha$  ELISA kit will enable determination of IFN- $\alpha$  levels in serum, plasma, and tissue culture media. As such, it should prove an important tool in virology, immunomodulation, and immunotoxicology studies conducted in pigs.

## Ordering Information

[Order Online »](#)

PBL-47100-1	96 wells
-------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Quantifies Porcine IFN-Alpha (Pig IFN- $\alpha$ ) in **serum, plasma**, and tissue culture media (TCM)
- Sensitive detection to as low as **2.34 pg/ml**
- High precision reproducibility & reliability with inter-assay CV  $\leq$  10% and intra-assay CV  $\leq$  8%

## 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 Pig IFN Alpha in serum, plasma, and tissue culture media

Assay Time Incubation time: 3 hours 15 minutes

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

Crossreactivity No cross-reactivity was observed when using this kit to detect the following, each at up to 20 ng/ml – Pig: IFN-β, IFN-γ; Human: IFN-α, IFN-β, IFN-γ, IFN-ω; Mouse: IFN-α, IFN-β, IFN-γ; Rat: IFN-α

Sensitivity 2.34 – 150 pg/ml

Species Reactivity Porcine



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)