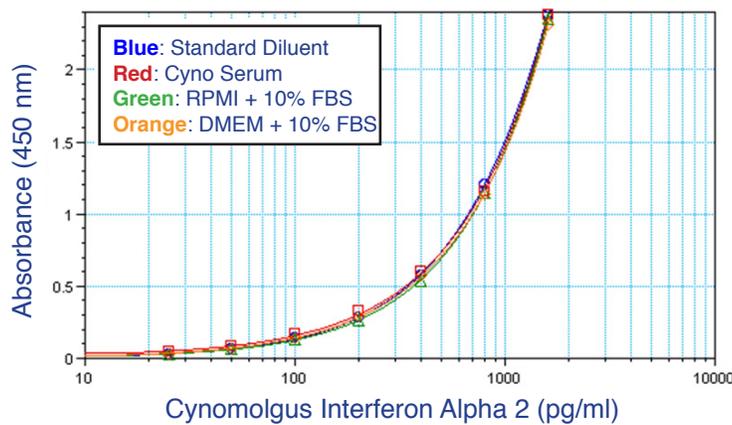




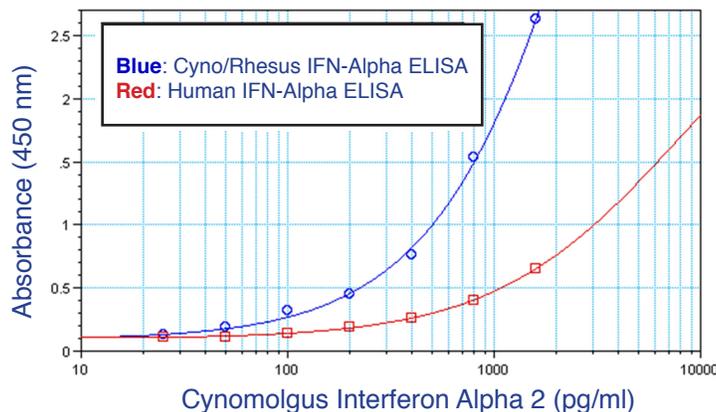
# VeriKine™ Cynomolgus/Rhesus Interferon Alpha Serum ELISA

Catalog No. 46100

- Enables sensitive detection of IFN- $\alpha$  to as low as 25 pg/ml
- Reproducible results with excellent inter- and intra-assay % CV (<15%)
- Quantifies IFN- $\alpha$ 2 in serum, plasma and tissue culture media
- Provides greater NHP IFN- $\alpha$  detection sensitivity and matrix compatibility than traditional human IFN- $\alpha$  ELISA kits
- Provides strong specificity for cynomolgus and rhesus IFN- $\alpha$ 2
- No cross-reactivity detected with human IFN- $\gamma$ , mouse IFN- $\alpha$  or rat IFN- $\alpha$



**Figure 1.** Standard Curve in Different Sample Matrices



**Figure 2.** Specifically designed for detecting NHP IFN- $\alpha$   
*Greater NHP IFN- $\alpha$  detection sensitivity and matrix compatibility*

Non-human primates (NHP) are considered a preferred model for the study of human diseases due to a close phylogenetic relationship to humans. The model is crucial to the elucidation of pathological mechanisms and may enable the development of novel diagnostic and therapeutic tools. Studies of innate and adaptive immunity, pharmacodynamics (PD), and immunotoxicology of interferon stimulating/inhibiting molecules, as well as Emerging Infectious Disease (EID) models, can all be further advanced with an interferon immunoassay that is designed specifically for NHP.

The interferon alpha gene family is comprised of multiple distinct genes that occupy a single chromosomal locus, and the congener proteins are produced and secreted by most somatic cells in response to pathogen exposure. Cynomolgus and Rhesus macaques have recently been shown to produce IFN alpha gene subtypes in response to Toll-like receptor agonist exposure.<sup>1</sup> The Cynomolgus/Rhesus IFN Alpha Serum ELISA has been developed to measure the presence of non-human primate IFN- $\alpha$ 2 in serum or tissue culture media samples by a sensitive sandwich enzyme linked immunosorbent assay (ELISA).

### Citations

1. Schramm, L., *et al.* (2012) High-Throughput Quantitative Real-Time Polymerase Chain Reaction Array for Absolute and Relative Quantification of Rhesus Macaque Types I, II and III Interferon and Their Subtypes. *Journal of Interferon & Cytokine Research* 32: 407-415.

### Product Information

Catalog Number	46100-1
Description	VeriKine™ Cynomolgus/Rhesus Interferon Alpha Serum ELISA
Size	1 x 96-well plate
Compatibility	Serum, Plasma, Cell Culture Supernatant
Assay Range	25 - 1600 pg/ml

For more information and product availability, please contact our customer service department by telephone (732-777-9123) or e-mail ([info@pbl assaysci.com](mailto:info@pbl assaysci.com)).



For more information, including a PBL white paper and product citations, visit our website.