



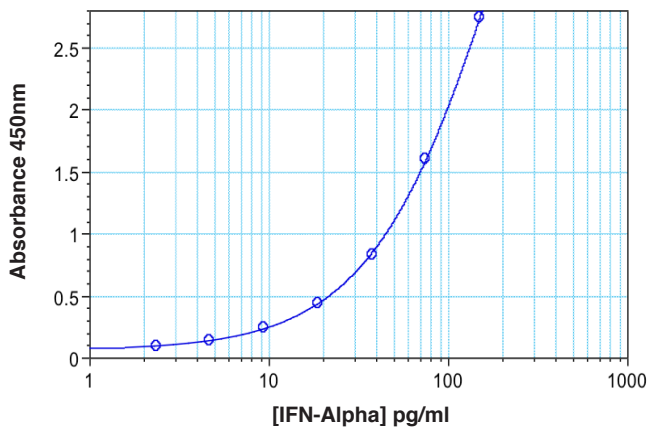
VeriKine-HS™

Pig Interferon Alpha ELISA

Catalog No. 47100

- Sensitive detection to as low as **2.34 pg/ml**
- Excellent matrix tolerance compatibility including **serum** and plasma
- Short incubation periods deliver accurate results in just over 3 hours
- High precision reproducibility with < 10% inter- and < 8% intra-assay CVs

Figure 1. Representative Standard Curve of Catalog No. 47100



The porcine model is gaining increasing interest from the scientific community as a non-rodent model for the study of human disease due to certain physiological similarities between swine and humans. This model may be used to study not only human diseases, but also viral and other diseases in livestock animals, especially as the transference of zoonotic viral diseases continues unabated. As infectious agents frequently cross species barriers and the probability of pandemic infections persists, better research tools are required.

This accurate and sensitive immunoassay enables detection of low levels of IFN-Alpha to advance porcine studies and help uncover insights into the pathologies of both human and livestock diseases, particularly in the area of novel viral infections which often induce interferon protein expression. This product could prove an important tool in virology, immunomodulation and immunotoxicology studies conducted in pigs.

pbl assay science

131 Ethel Road West, Suite 6 | Piscataway, NJ 08854

T (732) 777-9123 | **F** (732) 777-9141 | **E** info@pblassaysci.com | **W** www.pblassaysci.com

Product Information

Catalog Number	47100-1
Description	VeriKine-HS™ Pig Interferon Alpha ELISA
Size(s)	1 x 96-well plate
Compatibility	Serum, Plasma, Tissue Culture Media (TCM)
Assay Range	2.34 - 150 pg/ml

Visit our website for more information:

<https://www.pblassaysci.com/assay-kits/pig-ifn-alpha-elisa-kit-serum-plasma-tcm-47100>