

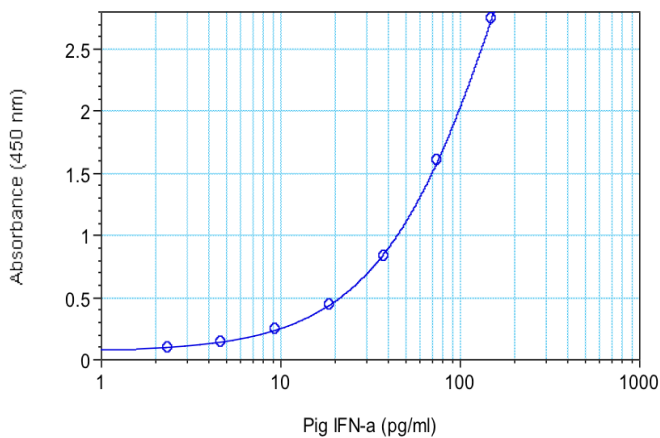


High Sensitivity Pig IFN Alpha ELISA

for sample in serum, plasma, and tissue culture media

Product Number: 47100-1

VeriKine-HS Pig IFN Alpha ELISA Kit



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

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.

PBL Assay Science is proud to announce the expansion of our porcine IFN product line to include a High Sensitivity Pig IFN Alpha ELISA kit. This accurate and sensitive assay enables detection of lower 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.