



## Product Notice

**Description:** VeriKine-HS™ Mouse Interferon Alpha All Subtype ELISA Kit

**Catalog No:** 42115-1

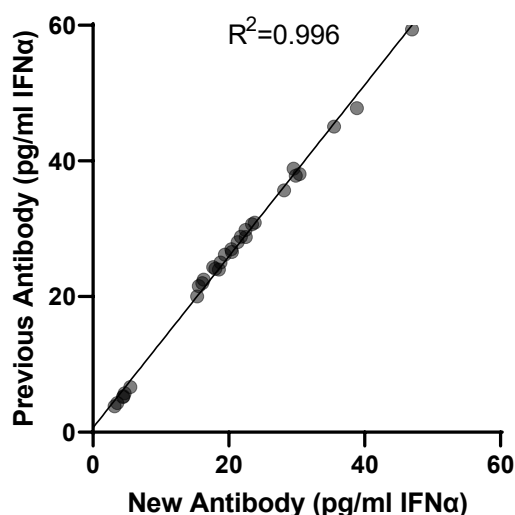
**Lots:** 7500 onward

**Date:** May 3, 2022

This notice is to inform you that a change has been made to the source of the detection antibody contained within the Antibody Concentrate component of the ELISA kit lots referenced above. The antibody stock used in the preparation of this component has been replaced with a new stock sourced from a different bleed from the same sheep donor. The affinity purification procedures were consistent between the antibody stock preparations, the formulation of the Antibody Concentrate component remains unchanged, and no other modifications have been made to the remaining kit components.

Please note that although the new Antibody Concentrate demonstrates a linear quantitative correlation with the previous one ( $R^2=0.996$ ), the new Antibody Concentrate yields lower relative sample concentrations. In a comparison study between antibody stocks, data generated in this kit with the provided Mouse IFN-Alpha 4 Standard has exhibited endogenous serum IFN-Alpha sample quantitation with the new antibody stock to be on average 77% of the concentration observed using the previous antibody. Please see the antibody correlation graph below:

**Correlation of Endogenous IFN $\alpha$  Measures  
using Two 42115 Antibody Concentrates (n=30)**



We thank you for your patience as PBL continues to improve the quality of its products. Please feel free to contact us at [techservice@pbl assaysci.com](mailto:techservice@pbl assaysci.com) if you have any product related questions.