

Product Notice

Description: VeriKine™ Cynomolgus/Rhesus Interferon Alpha Serum ELISA Kit

Catalog No: 46100-1

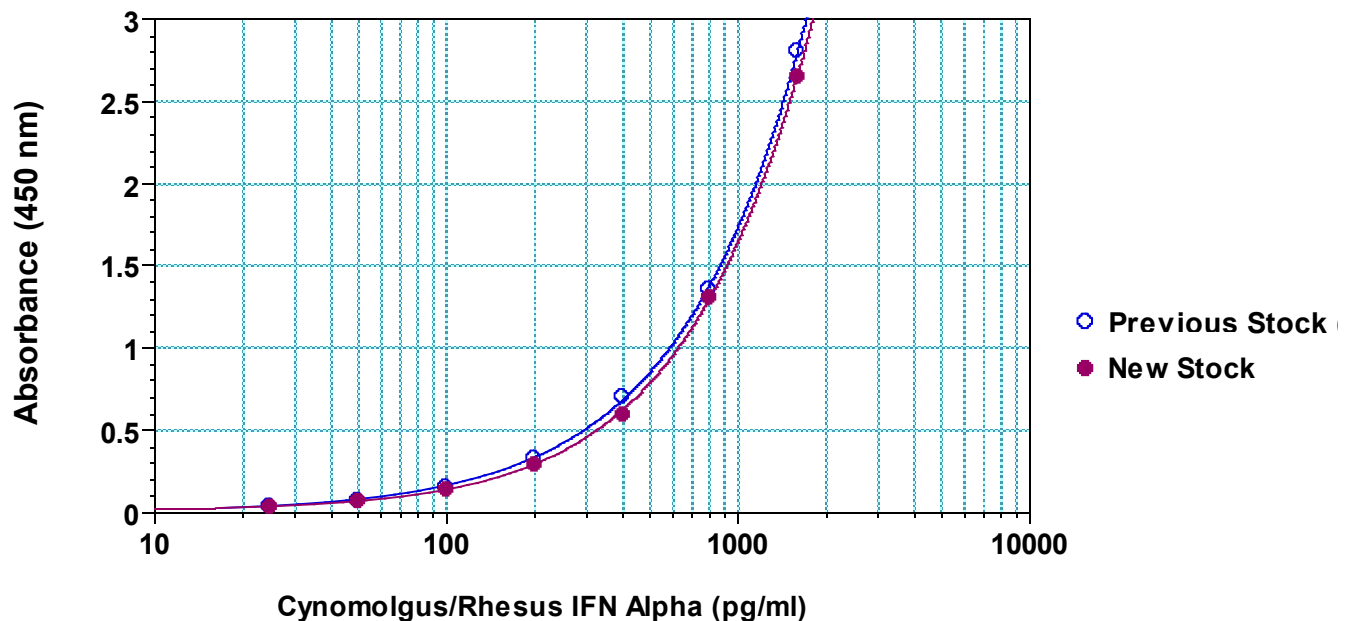
Lots: 7185 onward

Date: February 28, 2020

This notice is to inform you that a change has been made to the interferon standard used in the Cynomolgus/Rhesus Interferon Alpha Serum ELISA kit listed above. The formulation of the interferon standard remains the same, but the cynomolgus/rhesus interferon alpha stock used in preparation of this standard has been replaced with a new stock. No changes to the remaining kit components have been implemented.

Please find example data below from a standard study comparing the new cynomolgus/rhesus interferon alpha stock with the previous stock in the 46100 ELISA kit.

IFN α Standard Stock Comparison in 46100



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 if you have any product related questions.