

Product Notice

Description: VeriKine™ Human Interferon ELISA Kit

Catalog No: 41410-1, 41410-2

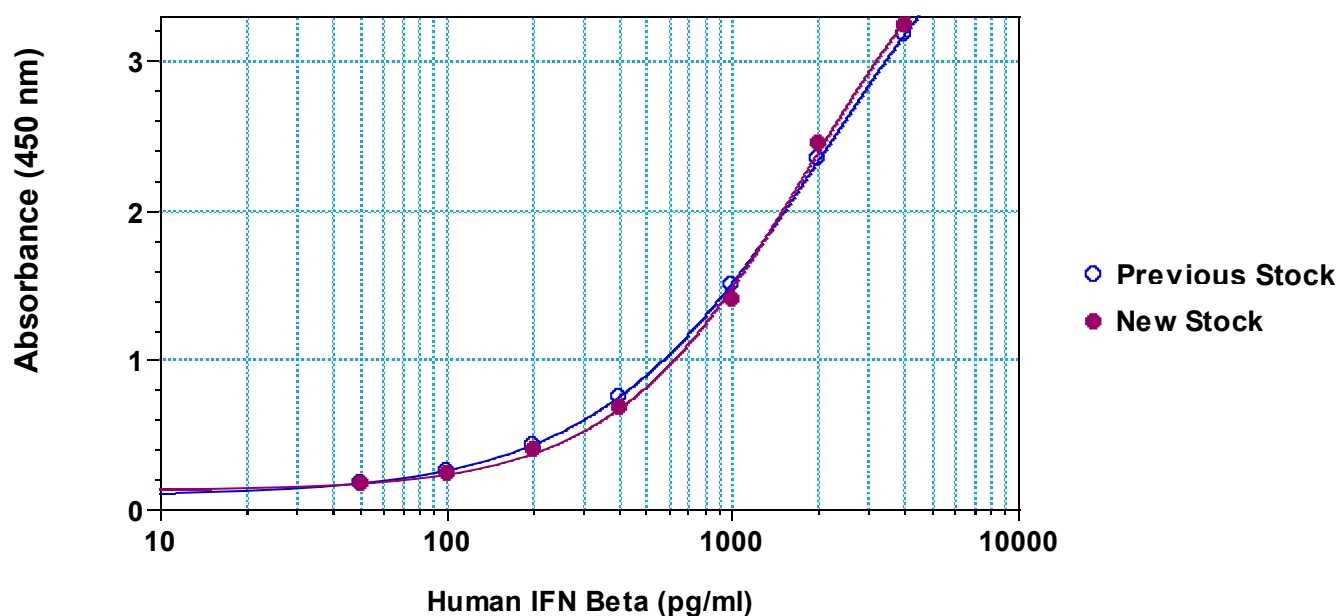
Lots: 6983 onward

Date: April 26, 2019

This notice is to inform you that a change has been made to the interferon standard used in the Human Interferon Beta ELISA kit listed above. The formulation of the interferon standard remains the same, but the human interferon beta 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 Human Interferon Beta stock with the previous stock in the 41410 ELISA kit.

IFN β Standard Stock Comparison in 41410



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.