Product Notice

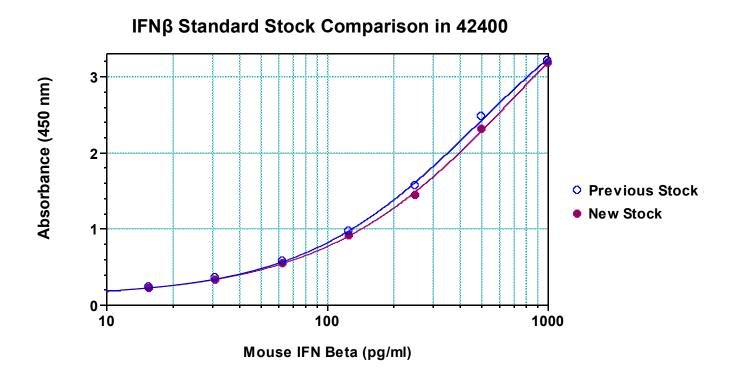
Description: VeriKine™ Mouse Interferon Beta ELISA Kit

Catalog No: 42400-1, 42400-2

Lots: 7047 onward **Date:** March 19, 2019

This notice is to inform you that a change has been made to the interferon standard used in the Mouse Interferon Beta ELISA kit listed above. The formulation of the interferon standard remains the same, but the mouse interferon beta stock used in preparation of this standard has been replaced with a new stock. The new standard has been shown to perform similarly to those derived from the previous source material. No changes to the remaining kit components have been implemented.

Please find data below from a standard study comparing the new Mouse Interferon Beta stock with the previous stock in the 42400 ELISA kit.



We thank you for your patience as PBL continues to improve the quality of its products. Please feel free to contact us at techservice@pblassaysci.com if you have any product related questions.