

Product Notice

Description: *Verikine*TM Mouse IFN Beta ELISA Kit

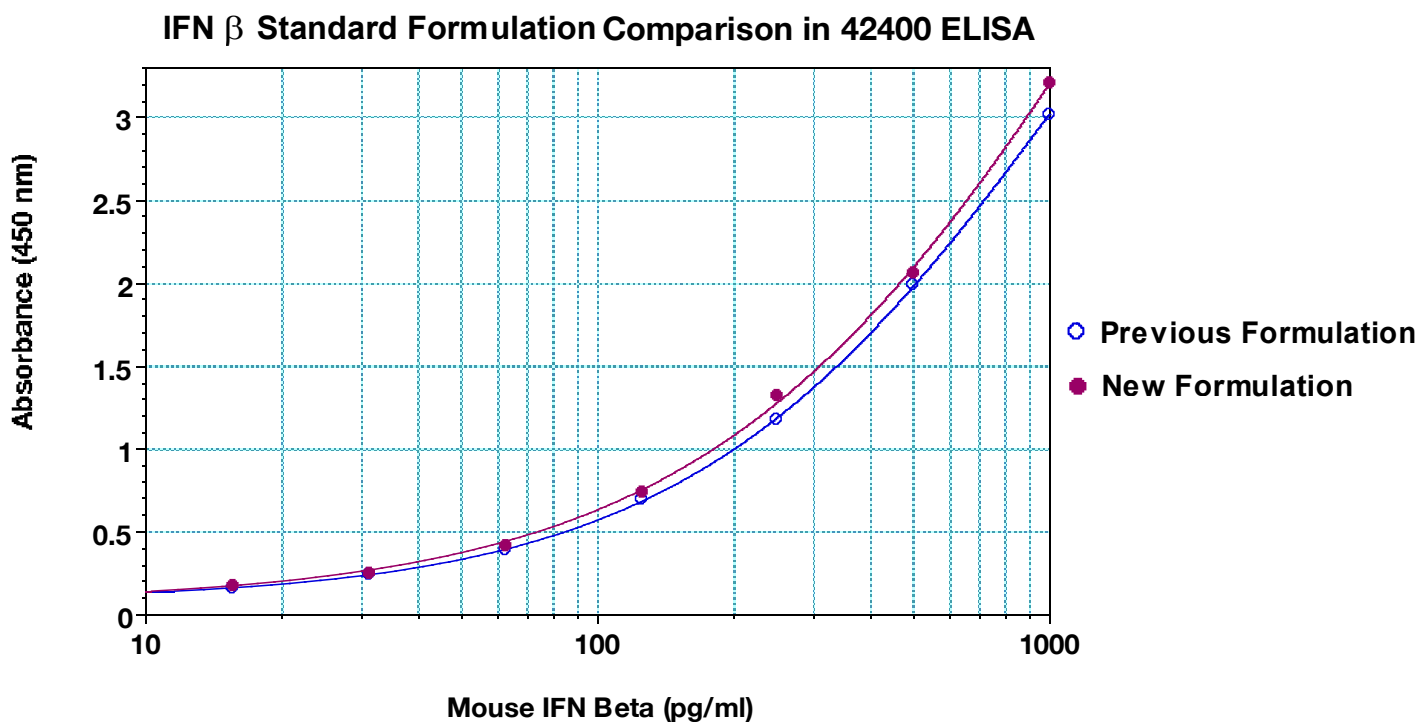
Catalog No: 42400-1, 42400-2

Lots: 6696 onward

Date: April 24, 2017

This notice is to inform you that there has been a formulation change to the above product. In order to improve stability and reduce variability, the buffer that the Mouse IFN Beta Standard component is stored in has been reformulated. This modification is intended to improve the long-term performance of the kit.

Please find data below from a study comparing the Mouse Interferon Beta Standard in the new buffer formulation with the previous buffer formulation 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@pbl assaysci.com if you have any product related questions.