



Product Notice

Date: 8/15/16

Catalog No: 41410-1, 41410-2

Lot: 6496 onward

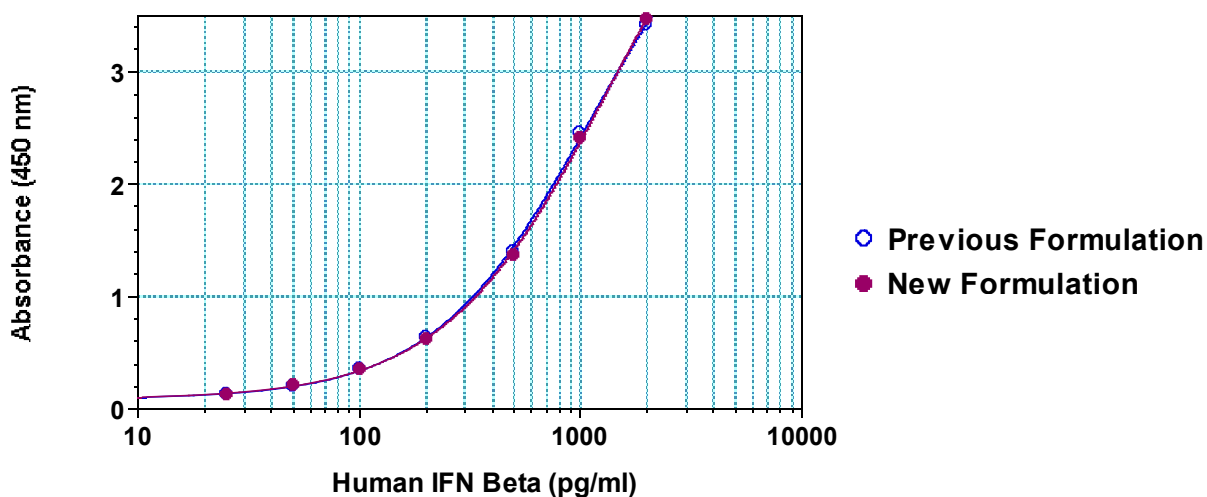
Description: Verikine Human Interferon Beta ELISA Kit

This notice is to inform you that there has been a formulation change to the Human Interferon Beta Standard in the above products. The Human Interferon Beta Standard component has been replaced with the standard formulation used in PBL's High Sensitivity Human Beta Serum ELISA kit (Cat. # 41415-1), designated with part number SMP146.

With this change, each individual 41410 kit will no longer be shipped in two separate containers . one containing the wet ice components and one containing the dry ice component (standard). **Beginning from lot 6496 onward, the entire 41410 kit, including the new standard formulation, will be shipped in a single container containing wet ice and can be stored at 4°C upon receipt.** As the new standard at 4°C will no longer be subject to the effect of freeze-thaw cycles, the standard will now be provided in a single vial suitable for multiple uses, rather than in three (prior 41410-1) or fifteen (prior 41410-2) separate, single-use vials. The concentration at which the new standard is provided remains at 100,000 pg/ml to match the prior standard and the serially diluted concentrations prepared in the standard curve also remain the same as in previous lots. Please refer to the protocol for modified instructions concerning handling of the protein during standard curve preparation and see the Certificate of Analysis for kit component information. All other components remain unchanged.

Please find data below from a study comparing the new Human Interferon Beta formulation SMP146 with the previous formulation SMP032 in the 41410 ELISA kit. The performance of each of these formulations has been shown to be highly similar in the 41410 ELISA:

IFN β Standard Formulation Comparison in 41410 ELISA



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.