

Product Notice

Date: 2/6/26

Catalog No: 41110-1, 41110-2

Lot: 7760, 7761, and onwards

Description: VeriKine™ Human Interferon Alpha Multi-Subtype Serum ELISA Kit

This notice is to inform you that a correction has been made to the instructions on the Certificate of Analysis & Protocol (CoA/Protocol) Rev.03 for the kits listed above.

The CoA/Protocol previously displayed some instructions calling for the standard or samples to be diluted in the Dilution Buffer component. However, the updated instructions direct the user to dilute the standard and samples in the relevant sample matrix or the Sample Diluent component. This change follows the original intention for the kit procedure reflected in earlier revisions.

Please be assured that these two diluents have similar formulations and results from side-by-side comparisons of standards and diluted samples prepared using both diluents have revealed insignificant differences in signal and quantification. This change should not affect the interpretation of results from assays prior to this notice.

All other information remains unchanged.

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.

Rev.00