



Product Notice

Description: VeriKine-HS™ Human Interferon Alpha All Subtype ELISA Kit, High Sensitivity (TCM)

Catalog No: 41135-1

Lots: 7501 onward

Date: September 8, 2022

This notice is to inform you that a modification has been made to the part number for the Antibody Concentrate component of the ELISA kit listed above. Originally listed as SMP322, the new part number that will be displayed on the bottle label and in the COA/Protocol will be SMP219.

Please note that this modification reflects a minor process change which better aligns internal procedures for improved efficiency. It does not indicate a formulation change or impact performance. No modifications have been made to the remaining kit components.

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

Rev. 00