



Certificate of Analysis

Protocol B

VeriKine-HS™ Human Interferon Beta Serum ELISA Kit
Certificate of Analysis for PROTOCOL B

Catalog No: 41415-1

Lot No: 7200

Expiration: April 30, 2021

Table with 4 columns: Kit Components, Part No., Lot No., Quantity. Lists components like Plate(s), Wash Solution Concentrate, and Assay Diluent.

Lot Specific Preparations

(Refer to VeriKine-HS™ Human Interferon Beta Serum ELISA Kit Protocol B for full instructions.)

Antibody Solution Preparation

Table for Antibody Solution Preparation with columns for Micro-plate Strips Used (2, 4, 6, 8, 10, 12) and rows for Antibody Concentrate and Assay Diluent.

HRP Solution Preparation

Table for HRP Solution Preparation with columns for Micro-plate Strips Used (2, 4, 6, 8, 10, 12) and rows for HRP Conjugate Concentrate and Assay Diluent.

Product Performance Specifications

Table for Product Performance Specifications with columns for Standard Diluent, Human Serum, and Tissue Culture Media (10% FBS) and rows for Intra-Assay CV and Inter-Assay CV.

Authorization

Released by: [Signature]

Date: May 13, 2020

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.