



Certificate of Analysis

Protocol A

VeriKine-HS™ Human Interferon Beta Serum ELISA Kit

Certificate of Analysis for **PROTOCOL A** (recommended protocol for improved serum performance)

Catalog No: 41415-1

Lot No: 7101

Expiration: July 15, 2020

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

Lot Specific Preparations

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

Antibody Solution Preparation

Table with 7 columns: Micro-plate Strips Used, and 6 columns for different strip counts (2, 4, 6, 8, 10, 12). Rows include Antibody Concentrate, Assay Diluent, Diluent Additive III, and Total Volume.

HRP Solution Preparation

Table with 7 columns: Micro-plate Strips Used, and 6 columns for different strip counts (2, 4, 6, 8, 10, 12). Rows include HRP Conjugate Concentrate and Assay Diluent.

Product Performance Specifications

Table with 4 columns: Standard Diluent, Human Serum, Tissue Culture Media (10% FBS). Rows include Intra-Assay CV and Inter-Assay CV.

Authorization

Released by: [Signature]

Date: August 13, 2019

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.