



### Certificate of Analysis

#### DIY Human IFN Lambda 3/1/2 (IL-28B/29/28A) ELISA Kit

Catalog No: 61840-1

Lot No: 7215

Expiration: March 15, 2024

Kit Components	Part No.	Lot No.	Quantity	Amount per vial
Capture Antibody	VDY066	KXU0419041	3	60 µg
Detection Antibody	VDY067	KLU0819041	3	24 µg
Standard	VDY068	1386478	3	120 ng
Streptavidin HRP	VDY803	P221868	3	2 ml

#### Lot Specific Solution Preparations

(Refer to **DIY Human IFN Lambda 3/1/2 (IL-28B/29/28A) ELISA Kit** Protocol for full instructions.)

Kit Components	Reagents Preparation	Working Concentration
Capture Antibody	Reconstitute with 0.5 ml of PBS	1.00 µg/ml
Detection Antibody	Reconstitute with 1.0 ml of Reagent Diluent	400 ng/ml
Standard	Reconstitute with 0.5 ml of Reagent Diluent	62.5-4000 pg/ml
Streptavidin HRP	Dilute with Reagent Diluent	40 -fold dilution

#### Authorization

Released by: \_\_\_\_\_

Date: April 14, 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.