



## Certificate of Analysis

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

**Catalog No:** 61830-1

**Lot No:**

**Expiration:**

Kit Components	Part No.	Lot No.	Quantity	Amount per vial
Capture Antibody	VDY896		1	µg
Detection Antibody	VDY897		1	µg
Standard	VDY898		3	ng
Streptavidin HRP	VDY803		1	ml

### Lot Specific Solution Preparations

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

Kit Components	Reagents Preparation	Working Concentration
Capture Antibody	Reconstitute with 1 ml of PBS	µg/ml
Detection Antibody	Reconstitute with 1.0 ml of Reagent Diluent	ng/ml
Standard	Reconstitute with 0.5 ml of Reagent Diluent	pg/ml
Streptavidin HRP	Dilute with Reagent Diluent	-fold dilution

### Authorization

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

Date:

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.

Rev. 00