

Certificate of Analysis

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

Catalog No: 61840-1

Lot No:

Expiration:

Kit Components	Part No.	Lot No.	Quantity	Amount per vial
Capture Antibody	VDY066		3	µg
Detection Antibody	VDY067		3	µg
Standard	VDY068		3	ng
Streptavidin HRP	VDY803		3	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	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. 01