



Certificate of Analysis

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

Catalog No: 61830-1

Lot No: 7242

Expiration: June 1, 2023

Kit Components	Part No.	Lot No.	Quantity	Amount per vial
Capture Antibody	VDY896	VJM0218051	1	360 µg
Detection Antibody	VDY897	KND0218051	1	18 µg
Standard	VDY898	1496198	3	145 ng
Streptavidin HRP	VDY803	P198628	1	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	2.00 µg/ml
Detection Antibody	Reconstitute with 1.0 ml of Reagent Diluent	100 ng/ml
Standard	Reconstitute with 0.5 ml of Reagent Diluent	125-8000 pg/ml
Streptavidin HRP	Dilute with Reagent Diluent	200 -fold dilution

Authorization

Released by: _____

Date: June 23, 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.