



Certificate of Analysis

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

Catalog No: 62830-1

Lot No: 7243

Expiration: June 1, 2023

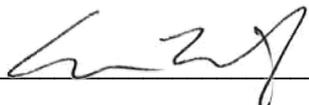
Kit Components	Part No.	Lot No.	Quantity	Amount per vial
Capture Antibody	VDY094	UCB0118111	3	60 µg
Detection Antibody	VDY095	UWZ0418111	3	15 µg
Standard	VDY096	1466490	6	40 ng
Streptavidin HRP	VDY803	P188503	3	2 ml

Lot Specific Solution Preparations

(Refer to **DIY Mouse IFN Lambda 2/3 (IL-28A/B) 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	250 ng/ml
Standard	Reconstitute with 0.5 ml of Reagent Diluent	31.2-2000 pg/ml
Streptavidin HRP	Dilute with Reagent Diluent	40 -fold dilution

Authorization

Released by: _____ 

Date: June 25, 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.