

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-15 Kit

Kit Lot Number: K0082331

Expiration Date: 31 Jul 2025

Kit Catalog Numbers: K151RD-Series; K156RD-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0047891	2-8°C	31 Dec 2025
Anti-hu IL-15 Antibody	D0081855	2-8°C	31 Oct 2025
Cytokine Panel 1 (human) Calibrator Blend	A0080331	2-8°C	31 Jul 2027
Diluent 3	M0030120	≤ -10°C	31 Jul 2025
Diluent 43	M0430053	≤ -10°C	31 Jan 2026

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision and Accuracy Specifications:

		Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-15	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-15	0.774	80% - 120%	< 20%	525	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-15	840	pg/mL

Dilution to Top of Curve = 1X

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		28 Nov 2023