

## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Kit Lot Number:** K0081894

**Expiration Date:** 29 Feb 2024

**Kit Catalog Numbers:** K151RD-Series; K156RD- Series

**Precision and Accuracy Specifications:**

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

Spot	Name	LLOQ			ULOQ		
		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	811	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0047616	2-8°C	30 Apr 2024
Anti-hu IL-15 Antibody	D0081610	2-8°C	31 Jul 2024
Cytokine Panel 1 (human) Calibrator Blend	A0080300	2-8°C	31 Jul 2026
Diluent 43	M0430046	≤ -10°C	29 Feb 2024
Diluent 3	M0030110	≤ -10°C	31 Mar 2024

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 FEB 2022