



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX PLUS Human IL-15 Kit
Lot Number: K0082713
Expiration Date: 30 SEP 2025
Catalog Number: K151RD-Series; K156RD-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-15 Antibody	D0082034	2-8°C	31 AUG 2026
Cytokine Panel 1 (human) Calibrator Blend	A0080300	2-8°C	31 JUL 2026
Human IL-15 Plate	Z0047851	2-8°C	30 SEP 2025
Diluent 3	M0030136	≤ -10°C	31 MAY 2027
Diluent 43	M0430056	≤ -10°C	30 SEP 2026

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

Precision, Accuracy, and Sensitivity 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

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 MAY 2025