

CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Kit Lot Number: K0082079

Expiration Date: 31 May 2024

Kit Catalog Numbers: K151RD-Series, K156RD-Series

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	811	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0047723	2-8°C	31 Dec 2024
Anti-hu IL-15 Antibody	D0081655	2-8°C	30 Sep 2024
Cytokine Panel 1 (human) Calibrator Blend	A0080300	2-8°C	31 Jul 2026
Diluent 3	M0030113	≤ -10°C	31 Jul 2024
Diluent 43	M0430047	≤ -10°C	31 May 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		24 AUG 2022