

**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-2 Kit

**Kit Lot Number:** K0082064

**Expiration Date:** 30 Jun 2024

**Kit Catalog Numbers:** K151QQ-Series; K156QQ-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-2	80% - 120%	< 20%	110% - 150%	< 20%	110% - 150%	< 20%	110% - 150%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-2	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	IL-2	1540	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-2 Plate	Z0047719	2-8°C	30 Nov 2024
Anti-hu IL-2 Antibody	D0081796	2-8°C	31 Aug 2025
Proinflammatory Panel 1 (human) Calibrator Blend	A0080328	2-8°C	30 Apr 2027
Diluent 2	M0020120	≤ -10°C	30 Jun 2024
Diluent 3	M0030114	≤ -10°C	30 Sep 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		21 SEP 2022