



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX PLUS Human IL-2 Kit
Lot Number: K0082459
Expiration Date: 30 JUN 2026
Catalog Number: K151QQ-Series; K156QQ-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-2 Antibody	D0082163	2-8°C	31 MAY 2027
Human IL-2 Plate	Z0047970	2-8°C	30 SEP 2026
Proinflammatory Panel 1 (human) Calibrator Blend	A0080385	2-8°C	31 DEC 2028
Diluent 2	M0020129	≤ -10°C	30 JUN 2026
Diluent 3	M0030127	≤ -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:

		Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-2	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-2	0.89	80% - 120%	< 20%	938	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-2	1380	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		30 MAY 2024