



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IL-2 Kit
Lot Number: K0082577
Expiration Date: 31 AUG 2026
Catalog Number: K151QQ-Series, K156QQ-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-2 Antibody	D0082004	2-8°C	31 AUG 2026
Human IL-2 Plate	Z0048035	2-8°C	31 MAR 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080411	2-8°C	31 OCT 2029
Diluent 2	M0020132	≤ -10°C	31 MAR 2027
Diluent 3	M0030132	≤ -10°C	31 MAR 2027

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

Precision, Accuracy, and Sensitivity 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%	100% - 150%	< 20%	100% - 150%	< 20%	100% - 150%	< 20%

Spot	Name	LLOQ			ULOQ		
		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	1410	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		10 DEC 2024