

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Human IL-2 Kit**Lot Number:** K0082893**Expiration Date:** 31 JAN 2028**Catalog Number:** K151QQ-Series, K156QQ-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-2 Antibody	D0082528	2-8°C	31 MAR 2029
Human IL-2, 96-Well Small Spot SECTOR Plate	Z0048230	2-8°C	31 JUL 2028
Proinflammatory Panel 1 (human) Calibrator Blend	A0080439	2-8°C	31 JUL 2030
Diluent 2	M0020136	≤ -10°C	31 JAN 2028
Diluent 3	M0030144	≤ -10°C	31 JAN 2028

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	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-2	1580	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		26 MAR 2026