



Extension Memo

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-2 Kit
Catalog Number(s): K151QQ-Series, K156QQ-Series
Lot Number: K0082610
New Expiration Date: 30 JUN 2025

Based on additional testing, the original date of expiration for this kit has been extended by 10 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-2 SECTOR Plate	Z0047671	31 AUG 2024	30 JUN 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		22 JAN 2025

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Human IL-2 Kit
Lot Number: K0082610
Expiration Date: 31 AUG 2024
Catalog Number: K151QQ-Series, K156QQ-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-2 Antibody	D0081796	2-8°C	31 AUG 2025
Human IL-2 SECTOR Plate	Z0047671	2-8°C	31 AUG 2024
Proinflammatory Panel 1 (human) Calibrator Blend	A0080400	2-8°C	30 JUN 2029
Diluent 2	M0020128	≤ -10°C	31 MAR 2026
Diluent 3	M0030129	≤ -10°C	30 NOV 2026

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	1430	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		20 JAN 2025