



Extension Memo


Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-2 Kit
Catalog Number(s): K152QQ-Series
Lot Number: K0082429
New Expiration Date: 30 Sep 2026

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Mouse IL-2 Plate	Z0047954	31 Jul 2026	30 Sep 2026
Diluent 41	M0410052	30 Sep 2025	30 Sep 2026
Diluent 45	M0450052	30 Sep 2025	30 Sep 2026
Anti-ms IL-2 Antibody	D0082002	31 Jul 2026	30 Sep 2026

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		31 Jul 2025

CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Kit Lot Number: K0082429

Expiration Date: 30 Sep 2025

Kit Catalog Numbers: K152QQ-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-2 Plate	Z0047954	2-8°C	31 Jul 2026
Anti-ms IL-2 Antibody	D0082002	2-8°C	31 Jul 2026
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080373	2-8°C	31 Aug 2028
Diluent 41	M0410052	≤ -10°C	30 Sep 2025
Diluent 45	M0450052	≤ -10°C	30 Sep 2025

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

Precision and Accuracy Specifications:

		Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-2	80% - 120%	< 20%	85% - 135%	< 25%	85% - 135%	< 25%	85% - 135%	< 25%


		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-2	1.03	75% - 125%	< 25%	1570	75% - 125%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-2	2580	pg/mL

Dilution to Top of Curve = 1X

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 Apr 2024