

**Extension Memo**

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-1 β Kit
Lot Number: K0082396
New Expiration Date: 30 NOV 2026
Catalog Number(s): K152QP-Series

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-1 β 96-Well Small Spot SECTOR Plate	Z0047930	30 APR 2026	30 NOV 2026
Anti-ms IL-1 β Antibody	D0081922	28 FEB 2026	30 NOV 2026
Diluent 41	M0410053	30 NOV 2025	30 NOV 2026
Diluent 45	M0450053	31 MAR 2026	30 NOV 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		24 Oct 2025

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-1 β Kit

Kit Lot Number: K0082396

Expiration Date: 30 Nov 2025

Kit Catalog Numbers: K152QP-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-1 β Plate	Z0047930	2-8°C	30 Apr 2026
Anti-ms IL-1 β Antibody	D0081922	2-8°C	28 Feb 2026
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080351	2-8°C	31 Dec 2027
Diluent 41	M0410053	$\leq$ -10°C	30 Nov 2025
Diluent 45	M0450053	$\leq$ -10°C	31 Mar 2026

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-1 β	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

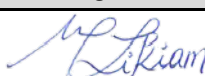
		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-1 β	0.720	75% - 125%	< 25%	1030	75% - 125%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-1 β	1740	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	Michelle Likiam	Quality		07 Dec 2023