



Extension Memo

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-1 β Kit
Catalog Number(s): K152QP-Series
Lot Number: K0082321
New Expiration Date: 31 Oct 2025

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
Anti-ms IL-1 β Antibody	D0081661	31 Oct 2024	31 Oct 2025
Proinflammatory Panel 1 (mouse) Calibrator	A0080252	30 Apr 2025	31 Oct 2025
Diluent 41	M0410052	30 Sep 2025	31 Oct 2025
Diluent 45	M0450052	30 Sep 2025	31 Oct 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		12 Dec 2024

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-1 β Kit
Kit Lot Number: K0082321
Expiration Date: 31 Oct 2024
Kit Catalog Numbers: K152QP-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-1 β	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		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	1530	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-1 β Plate	Z0047884	2-8°C	31 Dec 2025
Anti-ms IL-1 β Antibody	D0081661	2-8°C	31 Oct 2024
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080252	2-8°C	30 Apr 2025
Diluent 41	M0410052	\leq -10°C	30 Sep 2025
Diluent 45	M0450052	\leq -10°C	30 Sep 2025

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 AUG 2023