



### Extension Memo

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-1RA Kit  
**Catalog Number(s):** K151WT-Series  
**Lot Number:** K0082266  
**New Expiration Date:** 31 Jan 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
Human IL-1RA SECTOR Plate	Z0047844	30 Sep 2025	31 Jan 2026
Anti-hu IL-1RA Antibody	D0081720	31 Jan 2025	31 Jan 2026
Diluent 3	M0030120	31 Jul 2025	31 Jan 2026
Diluent 43	M0430050	31 Mar 2025	31 Jan 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		13 Dec 2024

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-1RA Kit  
**Kit Lot Number:** K0082266  
**Expiration Date:** 31 Jan 2025  
**Kit Catalog Numbers:** K151WT-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Cytokine Panel 2 (human) Control 1		Cytokine Panel 2 (human) Control 2		Cytokine Panel 2 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-1RA	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1RA	9.19	75% - 125%	< 25%	650	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	IL-1RA	1095	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1RA SECTOR Plate	Z0047844	2-8°C	30 Sep 2025
Anti-hu IL-1RA Antibody	D0081720	2-8°C	31 Jan 2025
Cytokine Panel 2 (human) Calibrator Blend	A0080295	2-8°C	31 May 2026
Diluent 3	M0030120	≤ -10°C	31 Jul 2025
Diluent 43	M0430050	≤ -10°C	31 Mar 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		02 JUN 2023