



## EXTENSION MEMO

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus Human IL-1 $\alpha$  Kit


**Catalog Number(s) / Size(s):** K151RB-Series

**Lot Number:** K0082101

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082101 has been extended to 28 FEB 2025.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-1 $\alpha$ Plate	Z0047717	30 Nov 2024	28 FEB 2025
Anti-hu IL-1 $\alpha$ Antibody	D0081571	30 Sep 2024	28 FEB 2025
Diluent 43	M0430046	29 Feb 2024	28 FEB 2025
Diluent 3	M0030110	31 Mar 2024	28 FEB 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		22 Feb 2024

This memo supersedes all other statements regarding date of expiration for this product.

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-1 $\alpha$  Kit

**Kit Lot Number:** K0082101

**Expiration Date:** 29 Feb 2024

**Kit Catalog Numbers:** K151RB-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-1 $\alpha$	80% - 120%	< 20%	95% - 135%	< 20%	95% - 135%	< 20%	95% - 135%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1 $\alpha$	2.85	80% - 120%	< 20%	278	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	IL-1 $\alpha$	448	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 $\alpha$ Plate	Z0047717	2-8°C	30 Nov 2024
Anti-hu IL-1 $\alpha$ Antibody	D0081571	2-8°C	30 Sep 2024
Cytokine Panel 1 (human) Calibrator Blend	A0080307	2-8°C	30 Sep 2026
Diluent 3	M0030110	$\leq$ -10°C	31 Mar 2024
Diluent 43	M0430046	$\leq$ -10°C	29 Feb 2024

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		06 SEP 2022