

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-1 $\alpha$  Kit  
**Kit Lot Number:** K0081600  
**Expiration Date:** 31 Aug 2022  
**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%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 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 $\alpha$	320	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 $\alpha$ Plate	Z0047426	2-8°C	30 Apr 2023
Anti-hu IL-1 $\alpha$ Antibody	D0081482	2-8°C	31 Dec 2023
Cytokine Panel 1 (human) Calibrator Blend	A0080240	2-8°C	30 Nov 2024
Diluent 43	M0430039	$\leq$ -10°C	30 Sep 2022
Diluent 3	M0030097	$\leq$ -10°C	31 Aug 2022

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		18 DEC 2020