

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-1RA Kit  
**Kit Lot Number:** K0081631  
**Expiration Date:** 31 Aug 2022  
**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	905	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1RA Plate	Z0047454	2-8°C	30 Jun 2023
Anti-hu IL-1RA Antibody	D0081241	2-8°C	31 Aug 2022
Cytokine Panel 2 (human) Calibrator Blend	A0080193	2-8°C	29 Feb 2024
Diluent 43	M0430040	≤ -10°C	30 Nov 2022
Diluent 3	M0030101	≤ -10°C	31 Dec 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		08 FEB 2021