

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-1RA Kit  
**Kit Catalog Numbers:** K151WT-Series  
**Kit Lot Number:** K0081265  
**Expiration Date:** 30 Sep 2021

**Acceptance Data/ Percent Recovery 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 Material Tested:**

Spot	Name	Item	Conc	Units
1	IL-1RA	Calibrator	905	pg/mL
		Cytokine Panel 2 (human) Control 1	514	pg/mL
		Cytokine Panel 2 (human) Control 2	93.1	pg/mL
		Cytokine Panel 2 (human) Control 3	54.4	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

IL-1RA - Spot 1		
Conc. pg/mL	ECL Signal	%CV
905	273,000	1.4
226	53,600	2.8
56.6	10,400	3.6
14.1	2,360	3.5
3.54	751	5.4
0.884	647	50.0
0.221	246	18.0
0.00	205	20.0

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1RA Plate	Z0047106	2-8°C	30 Nov 2021
Anti-hu IL-1RA Antibody	D0081093	2-8°C	31 Dec 2021
Cytokine Panel 2 (human) Calibrator Blend	A0080193	2-8°C	29 Feb 2024
Diluent 43	M0430032	≤ -10°C	30 Sep 2021
Diluent 3	M0030089	≤ -10°C	30 Sep 2021
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100 mL	Not Kit Specific
Cytokine Panel 2 (human) Control 1	Not Kit Specific
Cytokine Panel 2 (human) Control 2	Not Kit Specific
Cytokine Panel 2 (human) Control 3	Not Kit Specific

**Statement:**

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		07 AUG 2019