

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: K0081318
Expiration Date: 31 Dec 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	226,000	2.6
226	46,900	2.8
56.6	10,200	3.3
14.1	2,630	13.0
3.54	767	6.6
0.884	456	25.0
0.221	274	28.0
0.00	205	12.0

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1RA Plate	Z0047165	2-8°C	28 Feb 2022
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	M0430033	≤ -10°C	31 Jan 2022
Diluent 3	M0030090	≤ -10°C	31 Dec 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
Review/Approval	Karen Hamilla	Quality		02 OCT 2019