

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-1RA Kit
Kit Lot Number: K0081616
Expiration Date: 31 Dec 2021
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	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	M0430042	≤ -10°C	30 Apr 2023
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		04 JAN 2021