

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-1RA Kit
Kit Lot Number: K0082146
Expiration Date: 31 Oct 2024
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	1240	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1RA SECTOR Plate	Z0047774	2-8°C	31 Mar 2025
Anti-hu IL-1RA Antibody	D0081720	2-8°C	31 Jan 2025
Cytokine Panel 2 (human) Calibrator Blend	A0080251	2-8°C	30 Apr 2025
Diluent 3	M0030118	≤ -10°C	31 Mar 2025
Diluent 43	M0430049	≤ -10°C	31 Oct 2024

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		15 NOV 2022