

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-Plex Plus Human IL-8 Kit  
**Kit Lot Number:** K0081721  
**Expiration Date:** 31 May 2023  
**Kit Catalog Numbers:** K151RA-Series, K156RA-Series

**Precision and Accuracy Specifications:**

		Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-8	0.591	80% - 120%	< 20%	375	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	IL-8	602	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-8 Plate	Z0047527	2-8°C	30 Sep 2023
Anti-hu IL-8 Antibody	D0081399	2-8°C	30 Jun 2023
Proinflammatory Panel 1 (human) Calibrator Blend	A0080288	2-8°C	31 Mar 2026
Diluent 2	M0020115	≤ -10°C	31 May 2023
Diluent 3	M0030107	≤ -10°C	30 Sep 2023

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		02 JUL 2021