

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-8 Kit  
**Kit Lot Number:** K0082124  
**Expiration Date:** 30 Sep 2024  
**Kit Catalog Numbers:** K151RA-Series & K156RA-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-8	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		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	654	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-8 Plate	Z0047762	2-8°C	28 Feb 2025
Anti-hu IL-8 Antibody	D0081675	2-8°C	30 Sep 2024
Proinflammatory Panel 1 (human) Calibrator Blend	A0080339	2-8°C	31 Jul 2027
Diluent 2	M0020122	≤ -10°C	31 Oct 2024
Diluent 3	M0030116	≤ -10°C	30 Nov 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		21 OCT 2022