



### CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-8 Kit  
**Lot Number:** K0082544  
**Expiration Date:** 30 SEP 2026  
**Catalog Number:** K156RA-Series, K151RA-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-8 Antibody	D0082179	2-8°C	30 JUN 2027
Human IL-8 Plate	Z0048016	2-8°C	31 JAN 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080365	2-8°C	30 APR 2028
Diluent 2	M0020130	≤ -10°C	30 SEP 2026
Diluent 3	M0030128	≤ -10°C	31 OCT 2026

Additional components may be provided to execute the assay. Additional components are not lot-specific.

**Precision, Accuracy, and Sensitivity 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	593	pg/mL

Dilution to top of Curve: 1X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		30 SEP 2024