



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human IL-8 V2 Kit, SECTOR
Lot Number: K0082675
Expiration Date: 31 JUL 2027
Catalog Number: K151AWF-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-8 Antibody	D0082468	2-8°C	30 SEP 2028
Human IL-8 V2 96-Well Small Spot SECTOR Plate	Z0048099	2-8°C	30 SEP 2027
Diluent 58	M0580015	≤ -10°C	31 JUL 2027
Immune Response Panel 1 (human) Calibrator Blend	A0080422	≤ -70°C	30 SEP 2027

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

Precision, Accuracy, and Sensitivity Specifications:

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-8	0.716	80% - 120%	< 20%	585	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-8	4340	pg/mL

Dilution to top of Curve: 5X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 SEP 2025