



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IL-8 Kit
Lot Number: K0082997
Expiration Date: 30 APR 2028
Catalog Number: K151RA-Series, K156RA-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-8 Antibody	D0082360	2-8°C	30 APR 2028
Human IL-8, 96-Well Small Spot SECTOR Plate	Z0048288	2-8°C	31 DEC 2028
Proinflammatory Panel 1 (human) Calibrator Blend	A0080464	2-8°C	30 APR 2031
Diluent 2	M0020138	≤ -10°C	30 SEP 2028
Diluent 3	M0030148	≤ -10°C	31 JUL 2028

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

Precision, Accuracy, and Sensitivity 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%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 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	583	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		10 AUG 2026