



Extension Memo

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-8 Kit
Catalog Number(s): K151RA-Series; K156RA-Series
Lot Number: K0082419
New Expiration Date: 30 NOV 2025

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-8, 96-Well Small Spot SECTOR Plate	Z0047715	30 NOV 2024	30 NOV 2025
Diluent 2	M0020126	31 AUG 2025	30 NOV 2025
Diluent 3	M0030121	31 OCT 2025	30 NOV 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		23 JAN 2025



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IL-8 Kit
Lot Number: K0082419
Expiration Date: 30 NOV 2024
Catalog Number: K151RA-Series and K156RA-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-8 Antibody	D0081974	2-8°C	30 JUN 2026
Human IL-8, 96-Well Small Spot SECTOR Plate	Z0047715	2-8°C	30 NOV 2024
Proinflammatory Panel 1 (human) Calibrator Blend	A0080333	2-8°C	30 JUN 2027
Diluent 2	M0020126	≤ -10°C	31 AUG 2025
Diluent 3	M0030121	≤ -10°C	31 OCT 2025
MSD Read Buffer T (4X)	Y0140428	Room Temperature	30 JUN 2027

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	557	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		23 JAN 2025