



Extension Memo


Page 1 of 1

Product Description: V-PLEX® and V-PLEX PLUS Human IL-6 Kit
Catalog Number(s): K151QX-Series; K156QX-Series
Lot Number: K0082332
New Expiration Date: 28 Feb 2026

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-6 Plate	Z0047790	30 Apr 2025	28 Feb 2026
Anti-hu IL-6 Antibody	D0081822	31 Aug 2025	28 Feb 2026
Diluent 2	M0020124	28 Feb 2025	28 Feb 2026
Diluent 3	M0030118	31 Mar 2025	28 Feb 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		08 Apr 2025

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX PLUS Human IL-6 Kit
Kit Lot Number: K0082332
Expiration Date: 28 Feb 2025
Kit Catalog Numbers: K151QX-Series; K156QX-Series

Precision and Accuracy 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-6	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-6	0.633	80% - 120%	< 20%	488	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IL-6	725	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-6 Plate	Z0047790	2-8°C	30 Apr 2025
Anti-hu IL-6 Antibody	D0081822	2-8°C	31 Aug 2025
Proinflammatory Panel 1 (human) Calibrator Blend	A0080365	2-8°C	30 Apr 2028
Diluent 2	M0020124	≤ -10°C	28 Feb 2025
Diluent 3	M0030118	≤ -10°C	31 Mar 2025

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		27 JUL 2023