

**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-6 Kit

**Kit Lot Number:** K0082387

**Expiration Date:** 30 Nov 2025

**Kit Catalog Numbers:** K151QX-Series; K156QX-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-6 Plate	Z0047923	2-8°C	31 Mar 2026
Anti-hu IL-6 Antibody	D0082128	2-8°C	31 Jan 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080381	2-8°C	30 Sep 2028
Diluent 2	M0020127	≤ -10°C	30 Nov 2025
Diluent 3	M0030124	≤ -10°C	31 Mar 2026

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

**Precision and Accuracy 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-6	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%


Spot	Name	LLOQ			ULOQ		
		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	738	pg/mL

Dilution to Top of Curve = 1X

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		28 Nov 2023