

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX and V-PLEX Plus Human IL-6 Kit**Lot Number:** K0082925**Expiration Date:** 29 FEB 2028**Catalog Number:** K151QX-Series, K156QX-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-6 Antibody	D0082442	2-8°C	31 AUG 2028
Human IL-6, 96-Well Small Spot SECTOR Plate	Z0048249	2-8°C	30 SEP 2028
Proinflammatory Panel 1 (human) Calibrator Blend	A0080436	2-8°C	30 JUN 2030
Diluent 2	M0020137	≤ -10°C	29 FEB 2028
Diluent 3	M0030147	≤ -10°C	31 MAY 2028

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-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	733	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
<b>Review/Approval</b>	Karen Hamilla	Quality		06 MAY 2026