

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX Human IL-6 V2 Kit, SECTOR  
**Lot Number:** K0082648  
**Expiration Date:** 31 JUL 2027  
**Catalog Number:** K151AVE-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-6 Antibody	D0082463	2-8°C	31 AUG 2028
Human IL-6 V2 96-Well Small Spot SECTOR Plate	Z0048083	2-8°C	31 JUL 2027
Diluent 58	M0580015	≤ -10°C	31 JUL 2027
Cytolytic Panel 1 (human) Calibrator Blend	A0080419	≤ -70°C	31 JUL 2027

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

**Precision, Accuracy, and Sensitivity Specifications:**

Spot	Name	Calibrator		Cytolytic Panel 1 (human) Control 1		Cytolytic Panel 1 (human) Control 2		Cytolytic Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-6	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-6	0.436	80% - 120%	< 20%	780	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	IL-6	22100	pg/mL

Dilution to top of Curve: 20X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		22 AUG 2025