

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Mouse IL-6 Kit**Lot Number:** K0082771**Expiration Date:** 30 NOV 2026**Catalog Number:** K152QX-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms IL-6 Antibody	D0082248	2-8°C	30 SEP 2027
Mouse IL-6, 96-Well Small Spot SECTOR Plate	Z0048158	2-8°C	31 JAN 2028
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080401	2-8°C	30 JUN 2029
Diluent 41	M0410058	≤ -10°C	30 SEP 2027
Diluent 45	M0450054	≤ -10°C	30 NOV 2026

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 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-6	80% - 120%	< 20%	100% - 150%	< 25%	100% - 150%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-6	7.61	75% - 125%	< 25%	3140	75% - 125%	< 25%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	IL-6	4650	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	Michelle Likiam	Quality		10 OCT 2025