

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IL-6 Kit  
**Lot Number:** K0082835  
**Expiration Date:** 30 SEP 2027  
**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 Plate	Z0048208	2-8°C	31 MAY 2028
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080440	2-8°C	31 AUG 2030
Diluent 41	M0410059	≤ -10°C	31 DEC 2027
Diluent 45	M0450056	≤ -10°C	29 FEB 2028

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

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

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


		LLOQ			ULOQ		
Spot	Name	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	4950	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	Karen Hamilla	Quality		07 JAN 2026