



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IL-6 Kit  
**Lot Number:** K0082545  
**Expiration Date:** 28 FEB 2026  
**Catalog Number:** K152QX-Series

## Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-6 Antibody	D0081924	2-8°C	28 FEB 2026
Mouse IL-6 Plate	Z0048017	2-8°C	31 JAN 2027
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080351	2-8°C	31 DEC 2027
Diluent 41	M0410054	≤ -10°C	30 JUN 2026
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:

		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	4930	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		10 OCT 2024