

**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IL-6 Kit

**Kit Lot Number:** K0081621

**Expiration Date:** 31 Dec 2022

**Kit Catalog Numbers:** K152QX-Series

**Precision and Accuracy 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%	75% - 125%	< 25%	75% - 125%	< 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	5240	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-6 Plate	Z0047448	2-8°C	30 Jun 2023
Mouse IL-6 Antibody	D0081322	2-8°C	31 Dec 2022
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080264	2-8°C	31 Jul 2025
Diluent 41	M0410038	≤ -10°C	30 Jun 2023
Diluent 45	M0450040	≤ -10°C	31 Jul 2023

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		06 MAY 2021