

CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Kit Lot Number: K0081813

Expiration Date: 31 Jul 2023

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%

* See Additional Comments.

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	5070	pg/mL

Lot-Specific Components:


Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-6 Plate	Z0047569	2-8°C	31 Dec 2023
Anti-ms IL-6 Antibody	D0081443	2-8°C	31 Jul 2023
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080282	2-8°C	31 Mar 2026
Diluent 41	M0410041	≤ -10°C	30 Nov 2023
Diluent 45	M0450042	≤ -10°C	31 Dec 2023

Additional Comments:

Spot	Assay	Spot	Specification	Comment
1	IL-6	1	Control Conc.	Controls 1 & 2 are known to have biased % recovery (125-150%). Control 1 recovery value is 137.5%.

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		22 NOV 2021