

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Proinflammatory Panel 2 (rat) Kit  
**Kit Lot Number:** K0082346  
**Expiration Date:** 31 Oct 2024  
**Kit Catalog Numbers:** K15059-Series; K153QX-Series; K153QU-Series; K153OD-Series; K153QO-Series; K153QR-Series; K153QS-Series; K153QT-Series; K153QP-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Proinflammatory Panel 1 (rat) Control 1		Proinflammatory Panel 1 (rat) Control 2		Proinflammatory Panel 1 (rat) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- $\gamma$	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
2	IL-1 $\beta$	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
4	IL-4	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
5	IL-5	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
6	IL-6	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
7	KC/GRO	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
8	IL-10	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
9	IL-13	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
10	TNF- $\alpha$	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- $\gamma$	39.7	75% - 125%	< 25%	3750	75% - 125%	< 25%
2	IL-1 $\beta$	102	75% - 125%	< 25%	8100	75% - 125%	< 25%
4	IL-4	8.00	75% - 125%	< 25%	723	75% - 125%	< 25%
5	IL-5	82.0	75% - 125%	< 25%	3000	75% - 125%	< 25%
6	IL-6	96.9	75% - 125%	< 25%	8550	75% - 125%	< 25%
7	KC/GRO	21.7	75% - 125%	< 25%	728	75% - 125%	< 25%
8	IL-10	163	75% - 125%	< 25%	15700	75% - 125%	< 25%
9	IL-13	12.5	75% - 125%	< 25%	1080	75% - 125%	< 25%
10	TNF- $\alpha$	9.10	75% - 125%	< 25%	793	75% - 125%	< 25%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	IFN- $\gamma$	5287	pg/mL
2	IL-1 $\beta$	20400	pg/mL
4	IL-4	1133	pg/mL
5	IL-5	10153	pg/mL
6	IL-6	13900	pg/mL
7	KC/GRO	2683	pg/mL
8	IL-10	25900	pg/mL
9	IL-13	1693	pg/mL
10	TNF- $\alpha$	1283	pg/mL


**CERTIFICATE OF ANALYSIS**

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 2 (rat) Plate	Z0047824	2-8°C	31 Jul 2025
Anti-rat IFN- $\gamma$ Antibody	D0081993	2-8°C	31 Jul 2026
Anti-rat IL-1 $\beta$ Antibody	D0081995	2-8°C	31 Jul 2026
Anti-rat IL-4 Antibody	D0081994	2-8°C	31 Jul 2026
Anti-rat IL-5 Antibody	D0081996	2-8°C	31 Jul 2026
Anti-rat IL-6 Antibody	D0081997	2-8°C	31 Jul 2026
Anti-rat KC/GRO Antibody	D0081998	2-8°C	31 Jul 2026
Anti-rat IL-10 Antibody	D0081999	2-8°C	31 Jul 2026
Anti-rat IL-13 Antibody	D0082000	2-8°C	31 Jul 2026
Anti-rat TNF- $\alpha$ Antibody	D0082001	2-8°C	31 Jul 2026
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080376	2-8°C	31 Jul 2028
Blocker H	R0100022	Room Temperature	31 Oct 2024
Diluent 40	M0400019	$\leq -10^{\circ}\text{C}$	31 Oct 2024
Diluent 42	M0420018	$\leq -10^{\circ}\text{C}$	31 May 2025

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		12 SEP 2023