

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Proinflammatory Panel 1 (mouse) Kit  
**Kit Lot Number:** K0082254  
**Expiration Date:** 30 Jun 2025  
**Kit Catalog Numbers:** K15048-Series, K152QR-Series, K152QS-Series, K152QT-Series, K152QU-Series, K152QV-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	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%
3	IL-2	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-12p70	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- $\gamma$	0.390	75% - 125%	< 25%	570	75% - 125%	< 25%
2	IL-1 $\beta$	2.04	75% - 125%	< 25%	1030	75% - 125%	< 25%
3	IL-2	1.03	75% - 125%	< 25%	1570	75% - 125%	< 25%
4	IL-4	0.818	75% - 125%	< 25%	1060	75% - 125%	< 25%
5	IL-5	0.302	75% - 125%	< 25%	590	75% - 125%	< 25%
6	IL-6	7.61	75% - 125%	< 25%	3140	75% - 125%	< 25%
7	KC/GRO	3.29	75% - 125%	< 25%	1230	75% - 125%	< 25%
8	IL-10	7.26	75% - 125%	< 25%	2030	75% - 125%	< 25%
9	IL-12p70	179	75% - 125%	< 25%	20600	75% - 125%	< 25%
10	TNF- $\alpha$	0.980	75% - 125%	< 25%	403	75% - 125%	< 25%

**Non-Specific Binding Specifications:**

Spot	Name	NSB Spec
1	IFN- $\gamma$	< 1.0%
2	IL-1 $\beta$	< 1.0%
3	IL-2	< 1.0%
4	IL-4	< 1.0%
5	IL-5	< 1.0%
6	IL-6	< 1.0%
7	KC/GRO	< 1.0%
8	IL-10	< 1.0%
9	IL-12p70	< 1.0%
10	TNF- $\alpha$	< 1.0%

## CERTIFICATE OF ANALYSIS

### Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IFN- $\gamma$	1030	pg/mL
2	IL-1 $\beta$	1740	pg/mL
3	IL-2	2580	pg/mL
4	IL-4	1760	pg/mL
5	IL-5	1000	pg/mL
6	IL-6	4930	pg/mL
7	KC/GRO	2320	pg/mL
8	IL-10	3860	pg/mL
9	IL-12p70	27900	pg/mL
10	TNF- $\alpha$	622	pg/mL

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (mouse) Plate	Z0047836	2-8°C	31 Aug 2025
Anti-ms IFN- $\gamma$ Antibody	D0081921	2-8°C	28 Feb 2026
Anti-ms IL-1 $\beta$ Antibody	D0081922	2-8°C	28 Feb 2026
Anti-ms IL-2 Antibody	D0081828	2-8°C	31 Aug 2025
Anti-ms IL-4 Antibody	D0081829	2-8°C	31 Aug 2025
Anti-ms IL-5 Antibody	D0081923	2-8°C	31 Mar 2026
Anti-ms IL-6 Antibody	D0081924	2-8°C	28 Feb 2026
Anti-ms KC/GRO Antibody	D0081971	2-8°C	30 Jun 2026
Anti-ms IL-10 Antibody	D0081831	2-8°C	30 Jun 2026
Anti-ms IL-12p70 Antibody	D0081972	2-8°C	30 Jun 2026
Anti-ms TNF- $\alpha$ Antibody	D0081832	2-8°C	30 Nov 2025
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080351	2-8°C	31 Dec 2027
Diluent 41	M0410051	$\leq -10^{\circ}\text{C}$	31 Jul 2025
Diluent 45	M0450051	$\leq -10^{\circ}\text{C}$	30 Jun 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		18 May 2023