



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX PLUS Proinflammatory Panel 1 (mouse) Kit
Lot Number: K0082460
Expiration Date: 31 MAR 2026
Catalog Number: K15048-Series; K152QT-Series; K152QU-Series; K152QV-Series; K152QR-Series; K152QS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms IFN- γ Antibody	D0082064	2-8°C	31 OCT 2026
Anti-ms IL-10 Antibody	D0082067	2-8°C	31 OCT 2026
Anti-ms IL-12p70 Antibody	D0082068	2-8°C	31 OCT 2026
Anti-ms IL-1 β Antibody	D0082065	2-8°C	31 OCT 2026
Anti-ms IL-2 Antibody	D0082002	2-8°C	31 JUL 2026
Anti-ms IL-4 Antibody	D0082020	2-8°C	31 AUG 2026
Anti-ms IL-5 Antibody	D0082021	2-8°C	31 AUG 2026
Anti-ms IL-6 Antibody	D0082066	2-8°C	31 OCT 2026
Anti-ms KC/GRO Antibody	D0081971	2-8°C	30 JUN 2026
Anti-ms TNF- α Antibody	D0082022	2-8°C	31 AUG 2026
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080373	2-8°C	31 AUG 2028
Proinflammatory Panel 1 (mouse) Plate	Z0047971	2-8°C	30 SEP 2026
Diluent 41	M0410054	$\leq -10^{\circ}\text{C}$	30 JUN 2026
Diluent 45	M0450053	$\leq -10^{\circ}\text{C}$	31 MAR 2026

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

Precision, Accuracy, and Sensitivity 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- γ	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
2	IL-1 β	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- α	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

**CERTIFICATE OF ANALYSIS**

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	0.39	75% - 125%	< 25%	570	75% - 125%	< 25%
2	IL-1 β	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- α	0.98	75% - 125%	< 25%	403	75% - 125%	< 25%

Non-Specific Binding Specifications:


Spot	Name	Spec
1	IFN- γ	< 1.0%
2	IL-1 β	< 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- α	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	930	pg/mL
2	IL-1 β	1700	pg/mL
3	IL-2	2580	pg/mL
4	IL-4	1870	pg/mL
5	IL-5	923	pg/mL
6	IL-6	5160	pg/mL
7	KC/GRO	2150	pg/mL
8	IL-10	3770	pg/mL
9	IL-12p70	27000	pg/mL
10	TNF- α	610	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		19 JUN 2024