



**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX and V-PLEX Plus Proinflammatory Panel 1 (mouse) Kit  
**Lot Number:** K0082913  
**Expiration Date:** 30 NOV 2026  
**Catalog Number:** K15048-Series, K152QR-Series, K152QS-Series, K152QT-Series, K152QU-Series, K152QV-Series

**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms IFN- $\gamma$ Antibody	D0082443	2-8°C	31 AUG 2028
Anti-ms IL-10 Antibody	D0082511	2-8°C	31 JAN 2029
Anti-ms IL-1 $\beta$ Antibody	D0082444	2-8°C	31 AUG 2028
Anti-ms IL-4 Antibody	D0082510	2-8°C	28 FEB 2029
Anti-ms IL-6 Antibody	D0082448	2-8°C	31 AUG 2028
Anti-ms TNF- $\alpha$ Antibody	D0082452	2-8°C	30 NOV 2028
Mouse IL-12p70 Antibody	D0082508	2-8°C	28 FEB 2029
Mouse IL-2 Antibody	D0082506	2-8°C	28 FEB 2029
Mouse IL-5 Antibody	D0082507	2-8°C	28 FEB 2029
Mouse KC/GRO Antibody	D0082620	2-8°C	31 AUG 2029
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080440	2-8°C	31 AUG 2030
Proinflammatory Panel 1 (ms) 96-Well 10-Spot SECTOR Plate	Z0048162	2-8°C	31 JAN 2028
Diluent 41	M0410058	$\leq -10^{\circ}\text{C}$	30 SEP 2027
Diluent 45	M0450054	$\leq -10^{\circ}\text{C}$	30 NOV 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- $\gamma$	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%



### CERTIFICATE OF ANALYSIS

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	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%

#### Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- $\gamma$	991	pg/mL
2	IL-1 $\beta$	1640	pg/mL
3	IL-2	2610	pg/mL
4	IL-4	1710	pg/mL
5	IL-5	926	pg/mL
6	IL-6	4950	pg/mL
7	KC/GRO	2270	pg/mL
8	IL-10	3440	pg/mL
9	IL-12p70	31400	pg/mL
10	TNF- $\alpha$	602	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	Xiomara Talavera	Quality		23 MAR 2026