

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Cytokine Panel 1 (mouse) Kit  
**Kit Lot Number:** K0081727  
**Expiration Date:** 30 Apr 2023  
**Kit Catalog Numbers:** K15245-Series; K152XC-Series; K152XB-Series; K152WX-Series; K152QC-Series; K152RD-Series; K152WN-Series; K152NQ-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Cytokine Panel 1 (mouse) Control 1		Cytokine Panel 1 (mouse) Control 2		Cytokine Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-9	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	MCP-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	IL-33	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	IL-27p28/IL-30	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
5	IL-15	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	IL-17A/F	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
7	MIP-1 $\alpha$	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
8	IP-10	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
9	MIP-2	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-9	21.9	75% - 125%	< 20%	2600	75% - 125%	< 20%
2	MCP-1	4.42	75% - 125%	< 20%	325	75% - 125%	< 20%
3	IL-33	1.85	75% - 125%	< 20%	1950	75% - 125%	< 20%
4	IL-27p28/IL-30	5.91	75% - 125%	< 20%	6500	75% - 125%	< 20%
5	IL-15	43.2	75% - 125%	< 20%	26000	75% - 125%	< 20%
6	IL-17A/F	1.39	75% - 125%	< 20%	1620	75% - 125%	< 20%
7	MIP-1 $\alpha$	0.378	75% - 125%	< 20%	390	75% - 125%	< 20%
8	IP-10	6.58*	75% - 125%	< 20%	650	75% - 125%	< 20%
9	MIP-2	0.577	75% - 125%	< 20%	423	75% - 125%	< 20%

\*See Additional Comments

**Non-Specific Binding Specifications:**

Spot	Name	NSB Spec
1	IL-9	< 1.0%
2	MCP-1	< 1.0%
3	IL-33	< 1.0%
4	IL-27p28/IL-30	< 1.0%
5	IL-15	< 1.0%
6	IL-17A/F	< 1.0%
7	MIP-1 $\alpha$	< 1.0%
8	IP-10	< 1.0%
9	MIP-2	< 1.0%

**CERTIFICATE OF ANALYSIS**

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	IL-9	3645	pg/mL
2	MCP-1	496	pg/mL
3	IL-33	3025	pg/mL
4	IL-27p28/IL-30	9565	pg/mL
5	IL-15	39000	pg/mL
6	IL-17A/F	2805	pg/mL
7	MIP-1 $\alpha$	559	pg/mL
8	IP-10	1095	pg/mL
9	MIP-2	675	pg/mL

**Lot-Specific Components:**


Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (mouse) Plate	Z0047421	2-8°C	30 Apr 2023
Anti-ms IL-9 Antibody	D0081582	2-8°C	30 Jun 2024
Anti-ms MCP-1 Antibody	D0081583	2-8°C	30 Jun 2024
Anti-ms IL-33 Antibody	D0081584	2-8°C	30 Jun 2024
Anti-ms IL-27p28/IL-30 Antibody	D0081585	2-8°C	30 Jun 2024
Anti-ms IL-15 Antibody	D0081685	2-8°C	30 Sep 2024
Anti-ms IL-17A/F Antibody	D0081586	2-8°C	31 Jul 2024
Anti-ms MIP-1 $\alpha$ Antibody	D0081587	2-8°C	30 Jun 2024
Anti-ms IP-10 Antibody	D0081588	2-8°C	31 Jul 2024
Anti-ms MIP-2 Antibody	D0081589	2-8°C	31 Jul 2024
Cytokine Panel 1 (mouse) Calibrator Blend	A0080268	2-8°C	31 Aug 2025
Diluent 41	M0410038	$\leq -10^{\circ}\text{C}$	30 Jun 2023
Diluent 45	M0450040	$\leq -10^{\circ}\text{C}$	31 Jul 2023

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

**Additional Comments:**

Assay	Spot	Specification	Comment
IP-10	8	LLOQ Conc.	Did not meet specifications at 2.15 pg/mL; LLOQ reported at 6.58 pg/mL.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 JUN 2021