

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Gen. B Kit  
**Kit Lot Number:** K0082334  
**Expiration Date:** 31 Jul 2025  
**Kit Catalog Numbers:** K15705-series, K15712-series, K151ANG-series, K151ANQ-series, K151ANR-series, K156ANQ-series, K156ANR-series, K156ANG-series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Plate	Z0047893	2-8°C	31 Dec 2025
Human Eotaxin Antibody	D0082109	2-8°C	31 Jan 2027
Human MIP-1β Antibody	D0082110	2-8°C	28 Feb 2027
Human Eotaxin-3 Antibody	D0082111	2-8°C	31 Dec 2026
Human TARC Antibody	D0082112	2-8°C	31 Jan 2027
Human IP-10 Antibody	D0082113	2-8°C	31 Jan 2027
Human MIP-1α Antibody	D0082114	2-8°C	31 Jan 2027
Human MCP-1 Antibody	D0082115	2-8°C	31 Jan 2027
Human MDC Antibody	D0082116	2-8°C	31 Jan 2027
Human MCP-4 Antibody	D0082117	2-8°C	31 Jan 2027
Chemokine Panel 1 (hu) Gen. B Calibrator	A0080372	2-8°C	31 Jul 2028
Diluent 3	M0030123	≤ -10°C	31 Jan 2026
Diluent 57	M0570011	≤ -10°C	31 Jul 2025

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

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Chemokine Panel 1 (human) Gen. B Control 1		Chemokine Panel 1 (human) Gen. B Control 2		Chemokine Panel 1 (human) Gen. B Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Eotaxin	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	MIP-1β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	Eotaxin-3	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	TARC/CCL17	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	IP-10	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	MIP-1α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	MCP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	MDC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	MCP-4	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Eotaxin	2.14	80% - 120%	< 20%	910	80% - 120%	< 20%
2	MIP-1β	1.20	80% - 120%	< 20%	390	80% - 120%	< 20%
3	Eotaxin-3	11.1	80% - 120%	< 20%	11100	80% - 120%	< 20%
4	TARC/CCL17	2.14	80% - 120%	< 20%	481	75% - 125%	< 20%
5	IP-10	1.29	80% - 120%	< 20%	1240	80% - 120%	< 20%
6	MIP-1α	0.357	80% - 120%	< 20%	403	80% - 120%	< 20%
8	MCP-1	1.45	80% - 120%	< 20%	247	80% - 120%	< 20%
9	MDC	6.43	80% - 120%	< 20%	2930	75% - 125%	< 20%
10	MCP-4	0.518	80% - 120%	< 20%	234	80% - 120%	< 20%

## CERTIFICATE OF ANALYSIS

### Non-Specific Binding Specifications:


Spot	Name	NSB Spec
1	Eotaxin	< 1.0%
2	MIP-1 $\beta$	< 1.0%
3	Eotaxin-3	< 1.0%
4	TARC/CCL17	< 1.0%
5	IP-10	< 1.0%
6	MIP-1 $\alpha$	< 1.0%
8	MCP-1	< 1.0%
9	MDC	< 1.0%
10	MCP-4	< 1.0%

### Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Eotaxin	1450	pg/mL
2	MIP-1 $\beta$	594	pg/mL
3	Eotaxin-3	16900	pg/mL
4	TARC/CCL17	721	pg/mL
5	IP-10	1940	pg/mL
6	MIP-1 $\alpha$	618	pg/mL
8	MCP-1	397	pg/mL
9	MDC	4650	pg/mL
10	MCP-4	377	pg/mL

Dilution to Top of Curve = 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		02 FEB 2024