

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Kit

Kit Lot Number: K0082400

Expiration Date: 31 Jan 2026

Kit Catalog Numbers: K15047-Series; K15055-Series; K151NO-Series; K151NS-Series; K156NO-Series; K151NP-Series; K156NP-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0047934	2-8°C	31 May 2026
Anti-hu Eotaxin Antibody	D0081886	2-8°C	31 Jan 2026
Anti-hu MIP-1β Antibody	D0082160	2-8°C	31 May 2027
Anti-hu Eotaxin-3 Antibody	D0082077	2-8°C	30 Nov 2026
Anti-hu TARC Antibody	D0081888	2-8°C	31 Jan 2026
Anti-hu IP-10 Antibody	D0082078	2-8°C	30 Nov 2026
Anti-hu MIP-1α Antibody	D0081910	2-8°C	28 Feb 2026
Anti-hu IL-8 (HA) Antibody	D0081890	2-8°C	31 Jan 2026
Anti-hu MCP-1 Antibody	D0082159	2-8°C	31 May 2027
Anti-hu MDC Antibody	D0082079	2-8°C	30 Nov 2026
Human MCP-4 Antibody	D0081891	2-8°C	31 May 2026
Chemokine Panel 1 (human) Calibrator Blend	A0080369	2-8°C	30 Jun 2028
Diluent 3	M0030124	≤ -10°C	31 Mar 2026
Diluent 43	M0430055	≤ -10°C	30 Apr 2026

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) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) 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%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
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%
7	IL-8 (HA)	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	12.3	80% - 120%	< 20%	1120	80% - 120%	< 20%
2	MIP-1β	1.88	80% - 120%	< 20%	520	80% - 120%	< 20%
3	Eotaxin-3	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%
4	TARC/CCL17	3.32	80% - 120%	< 20%	1120	75% - 125%	< 20%
5	IP-10	1.37	80% - 120%	< 20%	500	80% - 120%	< 20%
6	MIP-1α	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%
7	IL-8 (HA)	713	75% - 125%	< 20%	43400	75% - 125%	< 20%
8	MCP-1	1.09	80% - 120%	< 20%	375	80% - 120%	< 20%
9	MDC	88.3	80% - 120%	< 20%	3700	75% - 125%	< 20%
10	MCP-4	1.49	80% - 120%	< 20%	472	80% - 120%	< 20%

CERTIFICATE OF ANALYSIS

Non-Specific Binding Specifications:


Spot	Name	NSB Spec
1	Eotaxin	< 1.0%
2	MIP-1 β	< 1.0%
3	Eotaxin-3	< 1.0%
4	TARC/CCL17	< 1.0%
5	IP-10	< 1.0%
6	MIP-1 α	< 1.0%
7	IL-8 (HA)	< 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	1860	pg/mL
2	MIP-1 β	1190	pg/mL
3	Eotaxin-3	6230	pg/mL
4	TARC/CCL17	1790	pg/mL
5	IP-10	2220	pg/mL
6	MIP-1 α	1170	pg/mL
7	IL-8 (HA)	74600	pg/mL
8	MCP-1	524	pg/mL
9	MDC	6060	pg/mL
10	MCP-4	718	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	Michelle Likiam	Quality		05 JAN 2024