



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX PLUS Chemokine Panel 1 (human) Kit
Lot Number: K0082447
Expiration Date: 30 APR 2026
Catalog Number: K15047-Series; K15055-Series; K151NO-Series; K156NO-Series; K156NP-Series; K151NP-Series; K151NS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu Eotaxin Antibody	D0082210	2-8°C	31 JUL 2027
Anti-hu Eotaxin-3 Antibody	D0082173	2-8°C	30 JUN 2027
Anti-hu IL-8 (HA) Antibody	D0082214	2-8°C	31 JUL 2027
Anti-hu IP-10 Antibody	D0082212	2-8°C	31 OCT 2027
Anti-hu MCP-1 Antibody	D0082215	2-8°C	31 AUG 2027
Anti-hu MDC Antibody	D0082175	2-8°C	30 JUN 2027
Anti-hu MIP-1 α Antibody	D0082036	2-8°C	31 AUG 2026
Anti-hu MIP-1 β Antibody	D0082211	2-8°C	31 JUL 2027
Anti-hu TARC Antibody	D0082174	2-8°C	30 JUN 2027
Chemokine Panel 1 (human) Calibrator Blend	A0080388	2-8°C	31 MAR 2029
Chemokine Panel 1 (human) Plate	Z0047965	2-8°C	30 SEP 2026
Human MCP-4 Antibody	D0081970	2-8°C	31 OCT 2026
Diluent 3	M0030128	$\leq -10^{\circ}\text{C}$	31 OCT 2026
Diluent 43	M0430055	$\leq -10^{\circ}\text{C}$	30 APR 2026

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

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Eotaxin	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	MIP-1 β	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-1 α	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%



CERTIFICATE OF ANALYSIS

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Eotaxin	12.3	80% - 120%	< 20%	1120	80% - 120%	< 20%
2	MIP-1 β	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-1 α	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%

Non-Specific Binding Specifications:

Spot	Name	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	< 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	1820	pg/mL
2	MIP-1 β	1170	pg/mL
3	Eotaxin-3	5790	pg/mL
4	TARC/CCL17	1750	pg/mL
5	IP-10	2320	pg/mL
6	MIP-1 α	1170	pg/mL
7	IL-8 (HA)	70900	pg/mL
8	MCP-1	553	pg/mL
9	MDC	5870	pg/mL
10	MCP-4	758	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		10 JUL 2024