

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Kit
Kit Lot Number: K0082149
Expiration Date: 31 Oct 2024
Kit Catalog Numbers: K15047-Series, K151NO-Series, K151NP-Series, K151NS-Series, K15055-Series, K156NP-Series, K156NO-Series

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	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	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	NSB Spec
1	Eotaxin	< 1.0%
2	MIP-1β	< 1.0%
3	Eotaxin-3	< 1.0%
4	TARC/CCL17	< 1.0%
5	IP-10	< 1.0%
6	MIP-1α	< 1.0%
7	IL-8	< 1.0%
8	MCP-1	< 1.0%
9	MDC	< 1.0%
10	MCP-4	< 1.0%

CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	Eotaxin	1770	pg/mL
2	MIP-1 β	1010	pg/mL
3	Eotaxin-3	6760	pg/mL
4	TARC/CCL17	1750	pg/mL
5	IP-10	2850	pg/mL
6	MIP-1 α	1180	pg/mL
7	IL-8	75400	pg/mL
8	MCP-1	564	pg/mL
9	MDC	5740	pg/mL
10	MCP-4	720	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0047776	2-8°C	31 Mar 2025
Anti-hu Eotaxin Antibody	D0081886	2-8°C	31 Jan 2026
Anti-hu MIP-1 β Antibody	D0081908	2-8°C	28 Feb 2026
Anti-hu Eotaxin-3 Antibody	D0081887	2-8°C	31 Jan 2026
Anti-hu TARC Antibody	D0081888	2-8°C	31 Jan 2026
Anti-hu IP-10 Antibody	D0081889	2-8°C	31 Jan 2026
Anti-hu MIP-1 α 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	D0081911	2-8°C	28 Feb 2026
Anti-hu MDC Antibody	D0081836	2-8°C	30 Sep 2025
Human MCP-4 Antibody	D0081891	2-8°C	31 May 2026
Chemokine Panel 1 (human) Calibrator Blend	A0080338	2-8°C	31 Aug 2027
Diluent 3	M0030118	\leq -10°C	31 Mar 2025
Diluent 43	M0430049	\leq -10°C	31 Oct 2024

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		06 JAN 2023