



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Kit**Lot Number:** K0082803**Expiration Date:** 28 FEB 2027**Catalog Number:** K15047-Series; K15055-Series; K156NO-Series; K151NO-Series; K151NP-Series; K151NS-Series; K156NP-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu Eotaxin Antibody	D0082472	2-8°C	30 SEP 2028
Anti-hu Eotaxin-3 Antibody	D0082474	2-8°C	30 SEP 2028
Anti-hu IL-8 (HA) Antibody	D0082478	2-8°C	31 OCT 2028
Anti-hu IP-10 Antibody	D0082476	2-8°C	31 OCT 2028
Anti-hu MCP-1 Antibody	D0082479	2-8°C	30 SEP 2028
Anti-hu MDC Antibody	D0082480	2-8°C	30 SEP 2028
Anti-hu MIP-1 α Antibody	D0082477	2-8°C	30 SEP 2028
Anti-hu MIP-1 β Antibody	D0082530	2-8°C	31 MAR 2029
Anti-hu TARC Antibody	D0082531	2-8°C	31 MAR 2029
Chemokine Panel 1 (human) Calibrator Blend	A0080432	2-8°C	30 APR 2030
Chemokine Panel 1 (human) Plate	Z0048182	2-8°C	31 MAR 2028
Human MCP-4 Antibody	D0082481	2-8°C	30 SEP 2028
Diluent 3	M0030141	$\leq -10^{\circ}\text{C}$	30 NOV 2027
Diluent 43	M0430058	$\leq -10^{\circ}\text{C}$	28 FEB 2027

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

Precision, Accuracy, and Sensitivity 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-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	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	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	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	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	1890	pg/mL
2	MIP-1 β	890	pg/mL
3	Eotaxin-3	6330	pg/mL
4	TARC	1780	pg/mL
5	IP-10	2520	pg/mL
6	MIP-1 α	1180	pg/mL
7	IL-8	76100	pg/mL
8	MCP-1	519	pg/mL
9	MDC	5600	pg/mL
10	MCP-4	746	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 NOV 2025