

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Gen. B Kit
Lot Number: K0082723
Expiration Date: 30 JUN 2027
Catalog Number: 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 (hu) Gen. B 96-Well 10-Spot SECTOR Plate	Z0048058	2-8°C	30 JUN 2027
Chemokine Panel 1 (hu) Gen. B Calibrator Blend	A0080431	2-8°C	31 MAR 2030
Human Eotaxin Antibody	D0082395	2-8°C	30 JUN 2028
Human Eotaxin-3 Antibody	D0082397	2-8°C	31 JUL 2028
Human IP-10 Antibody	D0082399	2-8°C	30 JUN 2028
Human MCP-1 Antibody	D0082401	2-8°C	30 JUN 2028
Human MCP-4 Antibody	D0082403	2-8°C	31 JUL 2028
Human MDC Antibody	D0082402	2-8°C	30 JUN 2028
Human MIP-1 α Antibody	D0082400	2-8°C	30 JUN 2028
Human MIP-1 β Antibody	D0082396	2-8°C	30 JUN 2028
Human TARC Antibody	D0082398	2-8°C	30 JUN 2028
Diluent 3	M0030139	$\leq -10^{\circ}\text{C}$	30 AUG 2027
Diluent 57	M0570020	$\leq -10^{\circ}\text{C}$	31 OCT 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) 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-1 β	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-1 α	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	2.14	80% - 120%	< 20%	910	80% - 120%	< 20%
2	MIP-1 β	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-1 α	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%

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%
8	MCP-1	< 1.0%
9	MDC	< 1.0%
10	MCP-4	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Eotaxin	1430	pg/mL
2	MIP-1 β	633	pg/mL
3	Eotaxin-3	17400	pg/mL
4	TARC/CCL17	730	pg/mL
5	IP-10	1980	pg/mL
6	MIP-1 α	590	pg/mL
8	MCP-1	370	pg/mL
9	MDC	4500	pg/mL
10	MCP-4	353	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		03 JUL 2025