

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Gen. B Kit, SECTOR
Lot Number: K0082842
Expiration Date: 30 SEP 2027
Catalog Number: K15705-Series; K15712-Series; K151ANG-Series; K151ANQ-Series; K151ANR-Series;
 K156ANQ-Series; K156ANR-Series; K156ANG-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (hu) Gen. B 96-Well 10-Spot SECTOR Plate	Z0048203	2-8°C	12 JUN 2028
Chemokine Panel 1 (human) 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	D0082557	2-8°C	31 MAY 2029
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 SEP 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

		LLOQ			ULOQ		
Spot	Name	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 FEB 2026