

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX PLUS Chemokine Panel 1 (human) Gen. B Kit, SECTOR
Lot Number: K0082970
Expiration Date: 30 JUN 2028
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 (human) Gen. B 96-Well 10-Spot SECTOR Plate	Z0048276	2-8°C	30 NOV 2028
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080462	2-8°C	31 MAR 2031
Human Eotaxin Antibody	D0082639	2-8°C	30 NOV 2029
Human Eotaxin-3 Antibody	D0082641	2-8°C	30 NOV 2029
Human IP-10 Antibody	D0082643	2-8°C	31 DEC 2029
Human MCP-1 Antibody	D0082645	2-8°C	30 NOV 2029
Human MCP-4 Antibody	D0082647	2-8°C	31 DEC 2029
Human MDC Antibody	D0082646	2-8°C	30 NOV 2029
Human MIP-1 α Antibody	D0082644	2-8°C	30 NOV 2029
Human MIP-1 β Antibody	D0082640	2-8°C	31 DEC 2029
Human TARC Antibody	D0082642	2-8°C	30 NOV 2029
Diluent 3	M0030148	$\leq -10^{\circ}\text{C}$	31 JUL 2028
Diluent 57	M0570022	$\leq -10^{\circ}\text{C}$	30 JUN 2028

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	1.2	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	1500	pg/mL
2	MIP-1 β	668	pg/mL
3	Eotaxin-3	17800	pg/mL
4	TARC/CCL17	784	pg/mL
5	IP-10	1970	pg/mL
6	MIP-1 α	683	pg/mL
8	MCP-1	413	pg/mL
9	MDC	4740	pg/mL
10	MCP-4	365	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		12 AUG 2026