

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX and V-PLEX Plus Chemokine Panel 1 (human) Kit**Lot Number:** K0082999**Expiration Date:** 30 JUN 2028**Catalog Number:** K15047-Series, K151NO-Series, K151NP-Series, K151NS-Series, K15055-Series, K156NO-Series, K156NP-Series, K151AS-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) 96-Well 10-Spot SECTOR Plate	Z0048284	2-8°C	31 DEC 2028
Chemokine Panel 1 (human) Calibrator Blend	A0080456	2-8°C	31 DEC 2030
Human Eotaxin Antibody	D0082656	2-8°C	31 JAN 2030
Human Eotaxin-3 Antibody	D0082658	2-8°C	31 DEC 2029
Human IL-8 (HA) Antibody	D0082662	2-8°C	31 DEC 2029
Human IP-10 Antibody	D0082660	2-8°C	31 DEC 2029
Human MCP-1 Antibody	D0082663	2-8°C	31 JAN 2030
Human MCP-4 Antibody	D0082665	2-8°C	31 DEC 2029
Human MDC Antibody	D0082664	2-8°C	31 JAN 2030
Human MIP-1 α Antibody	D0082661	2-8°C	31 JAN 2030
Human MIP-1 β Antibody	D0082657	2-8°C	31 JAN 2030
Human TARC Antibody	D0082659	2-8°C	31 DEC 2029
Diluent 3	M0030150	$\leq -10^{\circ}\text{C}$	31 OCT 2028
Diluent 43	M0430060	$\leq -10^{\circ}\text{C}$	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) 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/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%
7	IL-8 (HA)	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	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/CCL17	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 (HA)	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 (HA)	< 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	1860	pg/mL
2	MIP-1 β	889	pg/mL
3	Eotaxin-3	5620	pg/mL
4	TARC/CCL17	1780	pg/mL
5	IP-10	2310	pg/mL
6	MIP-1 α	1180	pg/mL
7	IL-8 (HA)	75200	pg/mL
8	MCP-1	580	pg/mL
9	MDC	5560	pg/mL
10	MCP-4	744	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		28 AUG 2026