



**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Gen. B Kit  
**Lot Number:** K0082578  
**Expiration Date:** 28 FEB 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	Z0048036	2-8°C	31 MAR 2027
Chemokine Panel 1 (hu) Gen. B Calibrator Blend	A0080389	2-8°C	30 APR 2029
Human Eotaxin Antibody	D0082326	2-8°C	29 FEB 2028
Human Eotaxin-3 Antibody	D0082328	2-8°C	31 MAR 2028
Human IP-10 Antibody	D0082330	2-8°C	29 FEB 2028
Human MCP-1 Antibody	D0082332	2-8°C	31 MAR 2028
Human MCP-4 Antibody	D0082334	2-8°C	31 MAR 2028
Human MDC Antibody	D0082333	2-8°C	29 FEB 2028
Human MIP-1 $\alpha$ Antibody	D0082359	2-8°C	30 APR 2028
Human MIP-1 $\beta$ Antibody	D0082327	2-8°C	29 FEB 2028
Human TARC Antibody	D0082358	2-8°C	30 APR 2028
Diluent 3	M0030135	$\leq -10^{\circ}\text{C}$	31 MAY 2027
Diluent 57	M0570017	$\leq -10^{\circ}\text{C}$	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) 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 $\beta$	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 $\alpha$	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 $\beta$	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 $\alpha$	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 $\beta$	< 1.0%
3	Eotaxin-3	< 1.0%
4	TARC/CCL17	< 1.0%
5	IP-10	< 1.0%
6	MIP-1 $\alpha$	< 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	1480	pg/mL
2	MIP-1 $\beta$	612	pg/mL
3	Eotaxin-3	15700	pg/mL
4	TARC/CCL17	778	pg/mL
5	IP-10	1800	pg/mL
6	MIP-1 $\alpha$	633	pg/mL
8	MCP-1	387	pg/mL
9	MDC	4710	pg/mL
10	MCP-4	374	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
<b>Review/Approval</b>	Karen Hamilla	Quality		24 JAN 2025