

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Kit
Kit Lot Number: K0081575
Expiration Date: 31 Jul 2022
Kit Catalog Numbers: K15047-Series, K15055-Series, K151NO-Series, K151NP-Series, K151NS-Series, K156NO-Series, K156NP-Series

Precision and Accuracy 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-1β	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	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	IP-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	MIP-1α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	IL-8	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%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Eotaxin	12.3	80% - 120%	< 20%	1120	80% - 120%	< 20%
2	MIP-1β	1.88	80% - 120%	< 20%	750	80% - 120%	< 20%
3	Eotaxin-3	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%
4	TARC	3.32	80% - 120%	< 20%	1120	80% - 120%	< 20%
5	IP-10	1.37	80% - 120%	< 20%	500	80% - 120%	< 20%
6	MIP-1α	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%
7	IL-8	713	75% - 125%	< 20%	13700	75% - 125%	< 20%
8	MCP-1	1.09	80% - 120%	< 20%	375	80% - 120%	< 20%
9	MDC	88.3	80% - 120%	< 20%	7500	80% - 120%	< 20%
10	MCP-4	10.0	75% - 125%	< 25%	469	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	NSB Spec
1	Eotaxin	< 1.0%
2	MIP-1β	< 1.0%
3	Eotaxin-3	< 1.0%
4	TARC	< 1.0%
5	IP-10	< 1.0%
6	MIP-1α	< 1.0%
7	IL-8	< 1.0%
8	MCP-1	< 1.0%
9	MDC	< 1.0%
10	MCP-4	< 1.0%

CERTIFICATE OF ANALYSIS

Page 2 of 2

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	Eotaxin	1870	pg/mL
2	MIP-1 β	1330	pg/mL
3	Eotaxin-3	5970	pg/mL
4	TARC	1830	pg/mL
5	IP-10	2360	pg/mL
6	MIP-1 α	1060	pg/mL
7	IL-8	64500	pg/mL
8	MCP-1	552	pg/mL
9	MDC	10200	pg/mL
10	MCP-4	724	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0047316	2-8°C	30 Sep 2022
Anti-hu Eotaxin Antibody	D0081287	2-8°C	31 Dec 2022
Anti-hu MIP-1 β Antibody	D0081181	2-8°C	31 Jul 2022
Anti-hu Eotaxin-3 Antibody	D0081260	2-8°C	31 Oct 2022
Anti-hu TARC Antibody	D0081261	2-8°C	31 Aug 2022
Anti-hu IP-10 Antibody	D0081262	2-8°C	31 Oct 2022
Anti-hu MIP-1 α Antibody	D0081508	2-8°C	30 Sep 2023
Anti-hu IL-8 (HA) Antibody	D0081264	2-8°C	31 Aug 2022
Anti-hu MCP-1 Antibody	D0081265	2-8°C	31 Aug 2022
Anti-hu MDC Antibody	D0081266	2-8°C	31 Aug 2022
Anti-hu MCP-4 Antibody	D0081267	2-8°C	31 Dec 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080228	2-8°C	31 Aug 2024
Diluent 43	M0430040	\leq -10°C	30 Nov 2022
Diluent 3	M0030101	\leq -10°C	31 Dec 2022

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality		16 NOV 2020