

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Cytokine Panel 1 (mouse) Kit
Kit Lot Number: K0081555
Expiration Date: 31 Mar 2022
Kit Catalog Numbers: K15245-Series, K152XC-Series, K152XB-Series, K152WX-Series, K152QC-Series, K152RD-Series, K152WN-Series, K152NQ-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (mouse) Control 1		Cytokine Panel 1 (mouse) Control 2		Cytokine Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-9	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	MCP-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	IL-33	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	IL-27p28/IL-30	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
5	IL-15	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	IL-17A/F	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
7	MIP-1 α	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
8	IP-10	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
9	MIP-2	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-9	21.9	75% - 125%	< 20%	2600	75% - 125%	< 20%
2	MCP-1	4.42	75% - 125%	< 20%	325	75% - 125%	< 20%
3	IL-33	1.85	75% - 125%	< 20%	1950	75% - 125%	< 20%
4	IL-27p28/IL-30	5.91	75% - 125%	< 20%	6500	75% - 125%	< 20%
5	IL-15	43.2	75% - 125%	< 20%	26000	75% - 125%	< 20%
6	IL-17A/F	1.39	75% - 125%	< 20%	1620	75% - 125%	< 20%
7	MIP-1 α	0.378	75% - 125%	< 20%	390	75% - 125%	< 20%
8	IP-10	2.15	75% - 125%	< 20%	650	75% - 125%	< 20%
9	MIP-2	0.577	75% - 125%	< 20%	423	75% - 125%	< 20%

Non-Specific Binding Specifications:

Spot	Name	NSB Spec
1	IL-9	< 1.0%
2	MCP-1	< 1.0%
3	IL-33	< 1.0%
4	IL-27p28/IL-30	< 1.0%
5	IL-15	< 1.0%
6	IL-17A/F	< 1.0%
7	MIP-1 α	< 1.0%
8	IP-10	< 1.0%
9	MIP-2	< 1.0%

CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IL-9	3720	pg/mL
2	MCP-1	447	pg/mL
3	IL-33	3010	pg/mL
4	IL-27p28/IL-30	8830	pg/mL
5	IL-15	39400	pg/mL
6	IL-17A/F	2380	pg/mL
7	MIP-1 α	739	pg/mL
8	IP-10	1100	pg/mL
9	MIP-2	559	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (mouse) Plate	Z0047357	2-8°C	30 Sep 2022
Anti-ms IL-9 Antibody	D0081497	2-8°C	31 Oct 2023
Anti-ms MCP-1 Antibody	D0081498	2-8°C	30 Nov 2023
Anti-ms IL-33 Antibody	D0081499	2-8°C	30 Nov 2023
Anti-ms IL-27p28/IL-30 Antibody	D0081500	2-8°C	30 Sep 2023
Anti-ms IL-15 Antibody	D0081297	2-8°C	30 Sep 2022
Anti-ms IL-17A/F Antibody	D0081502	2-8°C	31 Oct 2023
Anti-ms MIP-1 α Antibody	D0081503	2-8°C	30 Nov 2023
Anti-ms IP-10 Antibody	D0081504	2-8°C	30 Sep 2023
Anti-ms MIP-2 Antibody	D0081505	2-8°C	30 Nov 2023
Cytokine Panel 1 (mouse) Calibrator Blend	A0080218	2-8°C	31 Oct 2024
Diluent 41	M0410031	$\leq -10^{\circ}\text{C}$	30 Jun 2022
Diluent 45	M0450031	$\leq -10^{\circ}\text{C}$	31 Mar 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	Karen Hamilla	Quality		15 DEC 2020