



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Cytokine Panel 1 (mouse) Kit

Lot Number: K0082642

Expiration Date: 30 JUN 2026

Catalog Number: K15245-Series, K152XB-Series, K152XC-Series, K152NQ-Series, K152QC-Series, K152WX-Series, K152RD-Series, K152WN-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms IL-15 Antibody	D0082416	2-8°C	31 JUL 2028
Anti-ms IL-17A/F Antibody	D0082308	2-8°C	31 JAN 2028
Anti-ms IL-27p28/IL-30 Antibody	D0082306	2-8°C	31 JAN 2028
Anti-ms IL-33 Antibody	D0082305	2-8°C	31 MAR 2028
Anti-ms IL-9 Antibody	D0082303	2-8°C	31 JAN 2028
Anti-ms IP-10 Antibody	D0082310	2-8°C	31 MAR 2028
Anti-ms MCP-1 Antibody	D0082304	2-8°C	31 MAR 2028
Anti-ms MIP-1α Antibody	D0082309	2-8°C	31 JAN 2028
Anti-ms MIP-2 Antibody	D0082311	2-8°C	29 FEB 2028
Cytokine Panel 1 (mouse) 96-Well 10-Spot SECTOR Plate	Z0048010	2-8°C	31 JAN 2027
Cytokine Panel 1 (mouse) Calibrator Blend	A0080371	2-8°C	30 JUN 2028
Diluent 41	M0410054	≤ -10°C	30 JUN 2026
Diluent 45	M0450054	≤ -10°C	30 NOV 2026

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

Precision, Accuracy, and Sensitivity 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-1α	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-1α	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%



CERTIFICATE OF ANALYSIS

Non-Specific Binding Specifications:


Spot	Name	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%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-9	4160	pg/mL
2	MCP-1	574	pg/mL
3	IL-33	3220	pg/mL
4	IL-27p28/IL-30	9910	pg/mL
5	IL-15	41900	pg/mL
6	IL-17A/F	2890	pg/mL
7	MIP-1 α	630	pg/mL
8	IP-10	1030	pg/mL
9	MIP-2	692	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 FEB 2025