



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Cytokine Panel 1 (mouse) Kit**Lot Number:** K0082819**Expiration Date:** 30 NOV 2026**Catalog Number:** K15245-Series, K152RD-Series, K152WN-Series, K152WX-Series, K152XB-Series, K152XC-Series, K152NV-Series, K152NN-Series, K152NQ-Series, K152QC-Series, K152A0-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
Cytokine Panel 1 (mouse) Kit Calibrator	A0080371	2-8°C	30 JUN 2028
Cytokine Panel 1 (ms) 96-Well 10-Spot QuickPlex Ultra Plate	Z0048191	2-8°C	30 APR 2028
Mouse IL-15 Antibody	D0082457	2-8°C	31 OCT 2028
Mouse IL-17A/F Antibody	D0082458	2-8°C	30 SEP 2028
Mouse IL-27p28/IL-30 Antibody	D0082456	2-8°C	31 OCT 2028
Mouse IL-33 Antibody	D0082455	2-8°C	31 AUG 2028
Mouse IL-9 Antibody	D0082453	2-8°C	30 SEP 2028
Mouse IP-10 Antibody	D0082460	2-8°C	31 AUG 2028
Mouse MIP-1 α Antibody	D0082459	2-8°C	30 SEP 2028
Mouse MIP-2 Antibody	D0082461	2-8°C	30 NOV 2028
Mouse/Rat MCP-1 Antibody	D0082454	2-8°C	31 AUG 2028
Diluent 41	M0410056	$\leq -10^{\circ}\text{C}$	31 MAR 2027
Diluent 45	M0450054	$\leq -10^{\circ}\text{C}$	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-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%



CERTIFICATE OF ANALYSIS

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-9	21.9	75% - 125%	< 20%	2600	75% - 125%	< 20%
2	MCP-1	7.81	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%

*See Additional Comments

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



CERTIFICATE OF ANALYSIS

Additional Comments:

Spot	Name	Specification	Comment
8	IP-10	LLOQ CV	Specification for LLOQ is $\leq 20\%$, actual value 26.0% at target of 2.15pg/mL, was passing at 5.42pg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		29 JAN 2026