

**Extension Memo**

Product Description: V-PLEX® and V-PLEX PLUS Cytokine Panel 2 (human) Kit
Lot Number: K0082452
New Expiration Date: 31 MAR 2027
Catalog Number(s): K15084-Series; K151WP-Series; K151WO-Series; K151XC-Series;
K151WQ-Series; K151WZ Series; K151XE-Series

Based on additional testing, the original date of expiration for this kit has been extended by 11 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Cytokine Panel 2 (human) 96-Well 10-Spot	Z0047966	30 SEP 2026	31 MAR 2027
Diluent 3	M0030128	31 OCT 2026	31 MAR 2027
Diluent 43	M0430055	30 APR 2026	31 MAR 2027

Function	Name	Signature	Date
Quality	Karen Hamilla		19 Feb 2026



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX PLUS Cytokine Panel 2 (human) Kit
Lot Number: K0082452
Expiration Date: 30 APR 2026
Catalog Number: K15084-Series; K151WP-Series; K151WO-Series; K151XC-Series; K151WQ-Series; K151WZ-Series; K151XE-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-17A/F Antibody	D0082226	2-8°C	31 OCT 2027
Anti-hu IL-17B Antibody	D0082227	2-8°C	30 SEP 2027
Anti-hu IL-17C Antibody	D0082228	2-8°C	31 OCT 2027
Anti-hu IL-17D Antibody	D0082229	2-8°C	31 OCT 2027
Anti-hu IL-1RA Antibody	D0082230	2-8°C	30 SEP 2027
Anti-hu IL-3 Antibody	D0082231	2-8°C	30 SEP 2027
Anti-hu IL-9 Antibody	D0082232	2-8°C	30 SEP 2027
Anti-hu TSLP Antibody	D0082233	2-8°C	30 SEP 2027
Cytokine Panel 2 (human) Calibrator Blend	A0080346	2-8°C	31 DEC 2027
Cytokine Panel 2 (human) SECTOR Plate	Z0047966	2-8°C	30 SEP 2026
Diluent 3	M0030128	≤ -10°C	31 OCT 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026

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

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Cytokine Panel 2 (human) Control 1		Cytokine Panel 2 (human) Control 2		Cytokine Panel 2 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-17A/F	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	IL-17B	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	IL-17C	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	IL-17D	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	IL-1RA	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	IL-3	75% - 125%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	IL-9	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	TSLP	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-17A/F	7.57	80% - 120%	< 20%	3900	80% - 120%	< 20%
2	IL-17B	1.12	75% - 125%	< 20%	1040	75% - 125%	< 20%
3	IL-17C	3.84	75% - 125%	< 20%	1620	75% - 125%	< 20%
4	IL-17D	11.2	75% - 125%	< 20%	5200	75% - 125%	< 20%
6	IL-1RA	9.19	75% - 125%	< 25%	650	80% - 120%	< 20%
7	IL-3	12.6	75% - 125%	< 20%	1950	80% - 120%	< 20%
8	IL-9	2.23	80% - 120%	< 25%	975	80% - 120%	< 20%
10	TSLP	0.46	75% - 125%	< 20%	325	75% - 125%	< 20%



CERTIFICATE OF ANALYSIS

Non-Specific Binding Specifications:

Spot	Name	Spec
1	IL-17A/F	< 1.0%
2	IL-17B	< 1.0%
3	IL-17C	< 1.0%
4	IL-17D	< 1.0%
6	IL-1RA	< 1.0%
7	IL-3	< 1.0%
8	IL-9	< 1.0%
10	TSLP	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-17A/F	6000	pg/mL
2	IL-17B	1390	pg/mL
3	IL-17C	2510	pg/mL
4	IL-17D	8210	pg/mL
6	IL-1RA	1040	pg/mL
7	IL-3	3010	pg/mL
8	IL-9	1560	pg/mL
10	TSLP	475	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		21 JUN 2024