

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX and V-PLEX Plus Cytokine Panel 2 (human) Kit, SECTOR
Lot Number: K0082934
Expiration Date: 30 SEP 2027
Catalog Number: K151WN-Series, K151WO-Series, K151WP-Series, K151WQ-Series, K151WZ-Series, K151XC-Series, K151XE-Series, K15084-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
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) 96-Well 10-Spot SECTOR Plate	Z0048254	2-8°C	30 SEP 2028
Cytokine Panel 2 (human) Calibrator Blend	A0080444	2-8°C	30 SEP 2030
Diluent 3	M0030147	≤ -10°C	31 MAY 2028
Diluent 43	M0430059	≤ -10°C	30 SEP 2027

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 2 (human) Control 1		Cytokine Panel 2 (human) Control 2		Cytokine Panel 2 (human) Control 3	
		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%



CERTIFICATE OF ANALYSIS

Spot	Name	LLOQ			ULOQ		
		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%

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	6150	pg/mL
2	IL-17B	1230	pg/mL
3	IL-17C	2770	pg/mL
4	IL-17D	9010	pg/mL
6	IL-1RA	980	pg/mL
7	IL-3	3510	pg/mL
8	IL-9	1550	pg/mL
10	TSLP	521	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		11 JUN 2026