

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX and V-PLEX Plus Cytokine Panel 2 (human) Kit, QuickPlex**Lot Number:** K0082737**Expiration Date:** 30 NOV 2026**Catalog Number:** K15084-Series, K151WN-Series, K151WO-Series, K151WP-Series, K151WQ-Series, K151WT-Series, K151WZ-Series, K151XC-Series, K151XE-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-17A/F Antibody	D0082084	2-8°C	30 NOV 2026
Anti-hu IL-17B Antibody	D0082085	2-8°C	30 NOV 2026
Anti-hu IL-17C Antibody	D0082086	2-8°C	30 NOV 2026
Anti-hu IL-17D Antibody	D0082087	2-8°C	30 NOV 2026
Anti-hu IL-1RA Antibody	D0082088	2-8°C	31 JAN 2027
Anti-hu IL-3 Antibody	D0082089	2-8°C	30 NOV 2026
Anti-hu IL-9 Antibody	D0082500	2-8°C	31 DEC 2028
Anti-hu TSLP Antibody	D0082091	2-8°C	31 DEC 2026
Cytokine Panel 2 (hu) 96-Well 10-Spot QuickPlex Ultra Plate	Z0048118	2-8°C	30 SEP 2027
Cytokine Panel 2 (human) Calibrator Blend	A0080346	2-8°C	31 DEC 2027
Diluent 3	M0030140	≤ -10°C	30 SEP 2027
Diluent 43	M0430058	≤ -10°C	28 FEB 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%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 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%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
7	IL-3	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	IL-9	75% - 125%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	TSLP	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 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	80% - 120%	< 20%	1040	80% - 120%	< 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%	< 20%	650	75% - 125%	< 20%
7	IL-3	12.6	75% - 125%	< 25%	1950	80% - 120%	< 20%
8	IL-9	2.23	75% - 125%	< 20%	975	80% - 120%	< 20%
10	TSLP	0.46	80% - 120%	< 25%	325	80% - 120%	< 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	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		04 SEP 2025