

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX PLUS Cytokine Panel 2 (human) Kit
Kit Lot Number: K0082268
Expiration Date: 31 May 2024
Kit Catalog Numbers: K15084-Series; K151WN-Series; K151WO-Series; K151WP-Series; K151WQ-Series; K151WZ-Series; K151XC-Series; K151XE-Series

Precision and Accuracy 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%

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	0.460	75% - 125%	< 20%	325	75% - 125%	< 20%

Non-Specific Binding Specifications:

Spot	Name	NSB 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%

CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IL-17A/F	6500	pg/mL
2	IL-17B	1130	pg/mL
3	IL-17C	2645	pg/mL
4	IL-17D	8645	pg/mL
6	IL-1RA	1095	pg/mL
7	IL-3	3260	pg/mL
8	IL-9	1575	pg/mL
10	TSLP	502	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 2 (human) Plate	Z0047846	2-8°C	30 Sep 2025
Anti-hu IL-17A/F Antibody	D0082045	2-8°C	30 Sep 2026
Anti-hu IL-17B Antibody	D0081717	2-8°C	31 Jul 2025
Anti-hu IL-17C Antibody	D0081718	2-8°C	30 Jun 2025
Anti-hu IL-17D Antibody	D0081719	2-8°C	31 Mar 2025
Anti-hu IL-1RA Antibody	D0081961	2-8°C	31 May 2026
Anti-hu IL-3 Antibody	D0081559	2-8°C	31 May 2024
Anti-hu IL-9 Antibody	D0081722	2-8°C	31 Mar 2025
Anti-hu TSLP Antibody	D0081723	2-8°C	31 Jan 2025
Cytokine Panel 2 (human) Calibrator Blend	A0080295	2-8°C	31 May 2026
Diluent 3	M0030120	≤ -10°C	31 Jul 2025
Diluent 43	M0430050	≤ -10°C	31 Mar 2025

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 MAY 2023