

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Cytokine Panel 2 (human) Kit  
**Kit Lot Number:** K0081550  
**Expiration Date:** 30 Nov 2022  
**Kit Catalog Numbers:** K15084-Series, K151WO-Series, K151WP-Series, K151WQ-Series, K151WZ-Series, K151XC-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	6590	pg/mL
2	IL-17B	1790	pg/mL
3	IL-17C	2800	pg/mL
4	IL-17D	9050	pg/mL
6	IL-1RA	1240	pg/mL
7	IL-3	3710	pg/mL
8	IL-9	1620	pg/mL
10	TSLP	499	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 2 (human) Plate	Z0047394	2-8°C	28 Feb 2023
Anti-hu IL-17A/F Antibody	D0081554	2-8°C	31 Mar 2024
Anti-hu IL-17B Antibody	D0081555	2-8°C	31 Mar 2024
Anti-hu IL-17C Antibody	D0081556	2-8°C	31 May 2024
Anti-hu IL-17D Antibody	D0081557	2-8°C	30 Apr 2024
Anti-hu IL-1RA Antibody	D0081558	2-8°C	31 Mar 2024
Anti-hu IL-3 Antibody	D0081559	2-8°C	31 May 2024
Anti-hu IL-9 Antibody	D0081560	2-8°C	31 Mar 2024
Anti-hu TSLP Antibody	D0081561	2-8°C	31 Mar 2024
Cytokine Panel 2 (Human) Calibrator Blend	A0080251	2-8°C	30 Apr 2025
Diluent 43	M0430040	≤ -10°C	30 Nov 2022
Diluent 3	M0030101	≤ -10°C	31 Dec 2022

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		22 DEC 2020