



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Cytokine Panel 1 (human) Kit**Lot Number:** K0082806**Expiration Date:** 30 SEP 2027**Catalog Number:** K15050-Series, K151RC-Series, K151RE-Series, K15057-Series, K156RC-Series, K156RE-Series, K156RG-Series, K151RG-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu GM-CSF Antibody	D0082343	2-8°C	29 FEB 2028
Anti-hu IL-12/23p40 Antibody	D0082347	2-8°C	29 FEB 2028
Anti-hu IL-15 Antibody	D0082348	2-8°C	29 FEB 2028
Anti-hu IL-16 Antibody	D0082349	2-8°C	29 FEB 2028
Anti-hu IL-17A Antibody	D0082350	2-8°C	29 FEB 2028
Anti-hu IL-1 α Antibody	D0082344	2-8°C	29 FEB 2028
Anti-hu IL-5 Antibody	D0082345	2-8°C	29 FEB 2028
Anti-hu IL-7 Antibody	D0082346	2-8°C	29 FEB 2028
Anti-hu TNF- β Antibody	D0082351	2-8°C	29 FEB 2028
Anti-hu VEGF Antibody	D0082352	2-8°C	30 APR 2028
Cytokine Panel 1 (human) 96-Well 10-Spot SECTOR Plate	Z0048183	2-8°C	31 MAR 2028
Cytokine Panel 1 (human) Calibrator Blend	A0080435	2-8°C	31 MAY 2030
Diluent 3	M0030141	$\leq -10^{\circ}\text{C}$	30 NOV 2027
Diluent 43	M0430059	$\leq -10^{\circ}\text{C}$	30 SEP 2027

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

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	GM-CSF	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	IL-1 α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	IL-5	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	IL-7	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	IL-12/IL-23p40	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	IL-15	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	IL-16	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	IL-17A	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	TNF- β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	VEGF-A	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	GM-CSF	0.842	80% - 120%	< 20%	750	80% - 120%	< 20%
2	IL-1 α	2.85	80% - 120%	< 20%	278	80% - 120%	< 20%
3	IL-5	4.41	80% - 120%	< 20%	562	80% - 120%	< 20%
4	IL-7	0.851	80% - 120%	< 20%	563	80% - 120%	< 20%
5	IL-12/IL-23p40	1.32	80% - 120%	< 20%	2250	80% - 120%	< 20%
6	IL-15	0.774	80% - 120%	< 20%	525	80% - 120%	< 20%
7	IL-16	19.1	80% - 120%	< 20%	1870	80% - 120%	< 20%
8	IL-17A	3.19	80% - 120%	< 20%	3650	80% - 120%	< 20%
9	TNF- β	0.465	80% - 120%	< 20%	458	80% - 120%	< 20%
10	VEGF-A	7.70	80% - 120%	< 20%	562	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	Spec
1	GM-CSF	< 1.0%
2	IL-1 α	< 1.0%
3	IL-5	< 1.0%
4	IL-7	< 1.0%
5	IL-12/IL-23p40	< 1.0%
6	IL-15	< 1.0%
7	IL-16	< 1.0%
8	IL-17A	< 1.0%
9	TNF- β	< 1.0%
10	VEGF-A	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	GM-CSF	966	pg/mL
2	IL-1 α	435	pg/mL
3	IL-5	881	pg/mL
4	IL-7	878	pg/mL
5	IL-12/IL-23p40	3570	pg/mL
6	IL-15	889	pg/mL
7	IL-16	2560	pg/mL
8	IL-17A	5800	pg/mL
9	TNF- β	677	pg/mL
10	VEGF-A	1120	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		05 DEC 2025