

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Cytokine Panel 1 (human) Kit
Kit Lot Number: K0081628
Expiration Date: 30 Jun 2023
Kit Catalog Numbers: K15050-Series, K15057-Series, K151RC-Series, K151RG-Series, K156RG-Series, K156RC-Series, K151RE-Series, K156RE-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		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	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

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	0.546	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	7.70	80% - 120%	< 20%	562	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	NSB 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	< 1.0%

CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	GM-CSF	1280	pg/mL
2	IL-1 α	464	pg/mL
3	IL-5	906	pg/mL
4	IL-7	826	pg/mL
5	IL-12/IL-23p40	3650	pg/mL
6	IL-15	863	pg/mL
7	IL-16	3150	pg/mL
8	IL-17A	5340	pg/mL
9	TNF- β	705	pg/mL
10	VEGF	1520	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Plate	Z0047451	2-8°C	30 Jun 2023
Anti-hu GM-CSF Antibody	D0081570	2-8°C	31 Mar 2024
Anti-hu IL-1 α Antibody	D0081482	2-8°C	31 Dec 2023
Anti-hu IL-5 Antibody	D0081529	2-8°C	31 Dec 2023
Anti-hu IL-7 Antibody	D0081530	2-8°C	30 Jun 2024
Anti-hu IL-12/23p40 Antibody	D0081572	2-8°C	31 Mar 2024
Anti-hu IL-15 Antibody	D0081484	2-8°C	31 Oct 2023
Anti-hu IL-16 Antibody	D0081485	2-8°C	29 Feb 2024
Anti-hu IL-17A Antibody	D0081531	2-8°C	31 Dec 2023
Anti-hu TNF- β Antibody	D0081573	2-8°C	30 Jun 2024
Anti-hu VEGF Antibody	D0081574	2-8°C	31 Mar 2024
Cytokine Panel 1 (human) Calibrator Blend	A0080253	2-8°C	30 Apr 2025
Diluent 43	M0430043	$\leq -10^\circ\text{C}$	31 Jul 2023
Diluent 3	M0030104	$\leq -10^\circ\text{C}$	30 Jun 2023

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		05 MAY 2021