

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-Plex® and V-Plex Plus Proinflammatory Panel 1 (human) Kit  
**Kit Lot Number:** K0081744  
**Expiration Date:** 31 May 2023  
**Kit Catalog Numbers:** K151QR-Series; K151OD-Series; K15049-Series; K15052-Series; K15053-Series; K15056-Series; K15348-Series; K15346-Series; K15349-Series; K15345-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- $\gamma$	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	IL-1 $\beta$	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	IL-2	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	IL-4	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	IL-6	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	IL-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	IL-12p70	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	IL-13	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	TNF- $\alpha$	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- $\gamma$	1.76	80% - 120%	< 20%	938	80% - 120%	< 20%
2	IL-1 $\beta$	0.646	80% - 120%	< 20%	375	80% - 120%	< 20%
3	IL-2	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%
4	IL-4	0.218	80% - 120%	< 20%	158	80% - 120%	< 20%
5	IL-6	0.633	80% - 120%	< 20%	488	80% - 120%	< 20%
6	IL-8	0.591	80% - 120%	< 20%	375	80% - 120%	< 20%
7	IL-10	0.298	80% - 120%	< 20%	233	80% - 120%	< 20%
8	IL-12p70	1.22	80% - 120%	< 20%	315	80% - 120%	< 20%
9	IL-13	4.21	80% - 120%	< 20%	353	80% - 120%	< 20%
10	TNF- $\alpha$	0.690	80% - 120%	< 20%	248	80% - 120%	< 20%

**Non-Specific Binding Specifications:**

Spot	Name	NSB Spec
1	IFN- $\gamma$	< 1.0%
2	IL-1 $\beta$	< 1.0%
3	IL-2	< 1.0%
4	IL-4	< 1.0%
5	IL-6	< 1.0%
6	IL-8	< 1.0%
7	IL-10	< 1.0%
8	IL-12p70	< 1.0%
9	IL-13	< 1.0%
10	TNF- $\alpha$	< 1.0%

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IFN- $\gamma$	1530	pg/mL
2	IL-1 $\beta$	622	pg/mL
3	IL-2	1500	pg/mL
4	IL-4	225	pg/mL
5	IL-6	756	pg/mL
6	IL-8	617	pg/mL
7	IL-10	391	pg/mL
8	IL-12p70	464	pg/mL
9	IL-13	493	pg/mL
10	TNF- $\alpha$	337	pg/mL

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Plate	Z0047533	2-8°C	31 Oct 2023
Anti-hu IFN- $\gamma$ Antibody	D0081642	2-8°C	30 Sep 2024
Anti-hu IL-1 $\beta$ Antibody	D0081575	2-8°C	31 Aug 2024
Anti-hu IL-2 Antibody	D0081628	2-8°C	31 Aug 2024
Anti-hu IL-4 Antibody	D0081552	2-8°C	29 Feb 2024
Anti-hu IL-6 Antibody	D0081674	2-8°C	30 Sep 2024
Anti-hu IL-8 Antibody	D0081629	2-8°C	31 Aug 2024
Anti-hu IL-10 Antibody	D0081650	2-8°C	30 Sep 2024
Anti-hu IL-12p70 Antibody	D0081643	2-8°C	31 Aug 2024
Anti-hu IL-13 Antibody	D0081651	2-8°C	30 Sep 2024
Anti-hu TNF- $\alpha$ Antibody	D0081632	2-8°C	31 Jul 2024
Proinflammatory Panel 1 (human) Calibrator Blend	A0080289	2-8°C	31 Mar 2026
Diluent 2	M0020115	$\leq$ -10°C	31 May 2023
Diluent 3	M0030107	$\leq$ -10°C	30 Sep 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		01 JUL 2021