

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel1 (human) Kit
Kit Lot Number: K0082423
Expiration Date: 30 Jun 2026
Kit Catalog Numbers: K151QR-Series; K151OD-Series; K15049-Series; K15052-Series; K15053-Series; K15346-Series; K15349-Series; K15345-Series; K15348-Series; K15056-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Plate	Z0047950	2-8°C	31 Jul 2026
Anti-hu IFN- γ Antibody	D0082162	2-8°C	30 Apr 2027
Anti-hu IL-1 β Antibody	D0082107	2-8°C	31 Dec 2026
Anti-hu IL-2 Antibody	D0082163	2-8°C	31 May 2027
Anti-hu IL-4 Antibody	D0082063	2-8°C	30 Sep 2026
Anti-hu IL-6 Antibody	D0082128	2-8°C	31 Jan 2027
Anti-hu IL-8 Antibody	D0082108	2-8°C	31 Dec 2026
Anti-hu IL-10 Antibody	D0081975	2-8°C	30 Jun 2026
Anti-hu IL-12p70 Antibody	D0081984	2-8°C	31 Jul 2026
Anti-hu IL-13 Antibody	D0081976	2-8°C	30 Jun 2026
Anti-hu TNF- α Antibody	D0082164	2-8°C	30 Apr 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080382	2-8°C	30 Nov 2028
Diluent 2	M0020129	$\leq -10^\circ\text{C}$	30 Jun 2026
Diluent 3	M0030126	$\leq -10^\circ\text{C}$	31 Jul 2026

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

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- γ	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	IL-1 β	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- α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	1.76	80% - 120%	< 20%	938	80% - 120%	< 20%
2	IL-1 β	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- α	0.690	80% - 120%	< 20%	248	80% - 120%	< 20%

CERTIFICATE OF ANALYSIS

Non-Specific Binding Specifications:


Spot	Name	NSB Spec
1	IFN- γ	< 1.0%
2	IL-1 β	< 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- α	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	1340	pg/mL
2	IL-1 β	625	pg/mL
3	IL-2	1440	pg/mL
4	IL-4	218	pg/mL
5	IL-6	822	pg/mL
6	IL-8	586	pg/mL
7	IL-10	365	pg/mL
8	IL-12p70	478	pg/mL
9	IL-13	551	pg/mL
10	TNF- α	369	pg/mL

Dilution to Top of Curve = 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		21 MAR 2024