

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 1 (human) Kit
Kit Lot Number: K0082088
Expiration Date: 31 Oct 2024
Kit Catalog Numbers: K151OD-Series; K151QR-Series; K15049-Series; K15052-Series; K15053-Series; K15056-Series; K15345-Series; K15346-Series; K15348-Series; K15349-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- 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%

Non-Specific Binding Specifications:

Spot	Name	IFN- γ	IL-1 β	IL-2	IL-4	IL-6	IL-8	IL-10	IL-10	IL-12p70	IL-13	TNF- α
1	IFN- γ	NA	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%
2	IL-1 β	< 1.0%	NA	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%
3	IL-2	< 1.0%	< 1.0%	NA	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%
4	IL-4	< 1.0%	< 1.0%	< 1.0%	NA	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%
5	IL-6	< 1.0%	< 1.0%	< 1.0%	< 1.0%*	NA	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%
6	IL-8	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	NA	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%
7	IL-10	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	NA	NA	< 1.0%	< 1.0%	< 1.0%
8	IL-12p70	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	NA	< 1.0%	< 1.0%
9	IL-13	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	NA	< 1.0%
10	TNF- α	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	< 1.0%	NA

* See Additional Comments.

CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	1150	pg/mL
2	IL-1 β	613	pg/mL
3	IL-2	1540	pg/mL
4	IL-4	186	pg/mL
5	IL-6	730	pg/mL
6	IL-8	603	pg/mL
7	IL-10	370	pg/mL
8	IL-12p70	481	pg/mL
9	IL-13	491	pg/mL
10	TNF- α	375	pg/mL

Lot-Specific Components:


Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Plate	Z0047740	2-8°C	31 Jan 2025
Anti-hu IFN- γ Antibody	D0081808	2-8°C	30 Nov 2025
Anti-hu IL-1 β Antibody	D0081792	2-8°C	31 Aug 2025
Anti-hu IL-2 Antibody	D0081796	2-8°C	31 Aug 2025
Anti-hu IL-4 Antibody	D0081762	2-8°C	30 Apr 2025
Anti-hu IL-6 Antibody	D0081822	2-8°C	31 Aug 2025
Anti-hu IL-8 Antibody	D0081775	2-8°C	30 Jun 2025
Anti-hu IL-10 Antibody	D0081823	2-8°C	31 Aug 2025
Anti-hu IL-12p70 Antibody	D0081776	2-8°C	30 Jun 2025
Anti-hu IL-13 Antibody	D0081824	2-8°C	31 Oct 2025
Anti-hu TNF- α Antibody	D0081777	2-8°C	31 Aug 2025
Proinflammatory Panel 1 (human) Calibrator Blend	A0080328	2-8°C	30 Apr 2027
Diluent 2	M0020122	$\leq -10^{\circ}\text{C}$	31 Oct 2024
Diluent 3	M0030116	$\leq -10^{\circ}\text{C}$	30 Nov 2024

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

Additional Comments:

Assay	Spot	Specification	Comment
IL-6	5	NSB	MSD's upper limit for NSB for this panel is <1%; the actual measured NSB of Analyte IL-6 on the Analyte IL-4 spot was 1.046%

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

Name	Function	Signature	Date
Review/Approval Karen Hamilla	Quality		30 SEP 2022