



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX PLUS Proinflammatory Panel 1 (human) Kit
Lot Number: K0082511
Expiration Date: 31 MAR 2026
Catalog Number: 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
Anti-hu IFN- γ Antibody	D0082186	2-8°C	30 JUN 2027
Anti-hu IL-10 Antibody	D0082217	2-8°C	31 AUG 2027
Anti-hu IL-12p70 Antibody	D0082119	2-8°C	31 JAN 2027
Anti-hu IL-13 Antibody	D0082187	2-8°C	30 JUN 2027
Anti-hu IL-1 β Antibody	D0082176	2-8°C	30 JUN 2027
Anti-hu IL-2 Antibody	D0082234	2-8°C	31 OCT 2027
Anti-hu IL-4 Antibody	D0082177	2-8°C	30 JUN 2027
Anti-hu IL-6 Antibody	D0082249	2-8°C	30 SEP 2027
Anti-hu IL-8 Antibody	D0082179	2-8°C	30 JUN 2027
Anti-hu TNF- α Antibody	D0082218	2-8°C	31 AUG 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080400	2-8°C	30 JUN 2029
Proinflammatory Panel 1 (human) Plate	Z0048001	2-8°C	31 DEC 2026
Diluent 2	M0020128	$\leq -10^{\circ}\text{C}$	31 MAR 2026
Diluent 3	M0030129	$\leq -10^{\circ}\text{C}$	30 NOV 2026

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

Precision, Accuracy, and Sensitivity 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%

**CERTIFICATE OF ANALYSIS**

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	0.89	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	0.69	80% - 120%	< 20%	248	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	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	1060	pg/mL
2	IL-1 β	624	pg/mL
3	IL-2	1430	pg/mL
4	IL-4	183	pg/mL
5	IL-6	784	pg/mL
6	IL-8	575	pg/mL
7	IL-10	378	pg/mL
8	IL-12p70	442	pg/mL
9	IL-13	514	pg/mL
10	TNF- α	344	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		26 AUG 2024