

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Proinflammatory Panel 1 (rat) Kit**Lot Number:** K0082787**Expiration Date:** 28 FEB 2027**Catalog Number:** K15294-Series; K153J2-Series; K153K2-Series; K153L2-Series; K153M2-Series; K153N2-Series; K153P2-Series; K153Q2-Series; K153R2-Series; K153S2-Series; K153B2-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080410	2-8°C	30 JUN 2030
ProinflammPanel 1 (rt) 96-Well 10-Spot QuickPlex Ultra Plate	Z0048136	2-8°C	30 NOV 2027
Anti-rat IL-13 Antibody	D0082147	2-8°C	28 FEB 2027
Anti-rat TNF- α Antibody	D002497	2-8°C	30 NOV 2028
Rat IFN- γ Antibody	D0082488	2-8°C	30 NOV 2028
Rat IL-10 Antibody	D0082495	2-8°C	30 NOV 2028
Rat IL-1 β Antibody	D0082489	2-8°C	31 DEC 2028
Rat IL-2 Antibody	D0082490	2-8°C	31 DEC 2028
Rat IL-4 Antibody	D0082491	2-8°C	31 DEC 2028
Rat IL-5 Antibody	D0082492	2-8°C	31 DEC 2028
Rat IL-6 Antibody	D0082493	2-8°C	31 DEC 2028
Rat KC/GRO Antibody	D0082494	2-8°C	31 DEC 2028
Diluent 3	M0030140	$\leq -10^{\circ}\text{C}$	30 SEP 2027
Diluent 65	M0650003	$\leq -10^{\circ}\text{C}$	30 NOV 2027

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 (rat) Control 1		Proinflammatory Panel 1 (rat) Control 2		Proinflammatory Panel 1 (rat) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- γ	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	IL-1 β	80% - 120%	< 20%	70% - 130%	< 20%	70% - 130%	< 20%	70% - 130%	< 20%
3	IL-2	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	IL-4	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
5	IL-5	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	IL-6	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
7	KC/GRO	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
8	IL-10	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
9	IL-13	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
10	TNF- α	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%



CERTIFICATE OF ANALYSIS

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	1.21	75% - 125%	< 20%	292	75% - 125%	< 20%
2	IL-1 β	2.66	75% - 125%	< 20%	1230	75% - 125%	< 20%
3	IL-2	1.07	75% - 125%	< 20%	455	75% - 125%	< 20%
4	IL-4	3.62	75% - 125%	< 20%	357	75% - 125%	< 20%
5	IL-5	1.86	75% - 125%	< 20%	305	75% - 125%	< 20%
6	IL-6	188	75% - 125%	< 20%	40000	75% - 125%	< 20%
7	KC/GRO	2.23	75% - 125%	< 20%	604	75% - 125%	< 20%
8	IL-10	43.1	75% - 125%	< 25%	3900	75% - 125%	< 25%
9	IL-13	11.8	75% - 125%	< 25%	910	75% - 125%	< 25%
10	TNF- α	2.44	75% - 125%	< 25%	715	75% - 125%	< 25%

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-5	< 1.0%
6	IL-6	< 1.0%
7	KC/GRO	< 1.0%
8	IL-10	< 1.0%
9	IL-13	< 1.0%
10	TNF- α	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	438	pg/mL
2	IL-1 β	1530	pg/mL
3	IL-2	761	pg/mL
4	IL-4	535	pg/mL
5	IL-5	453	pg/mL
6	IL-6	94500	pg/mL
7	KC/GRO	997	pg/mL
8	IL-10	7150	pg/mL
9	IL-13	1460	pg/mL
10	TNF- α	1210	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	Michelle Likiam	Quality		09 OCT 2025