

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Proinflammatory Panel 1 (rat) Kit, SECTOR**Lot Number:** K0082558**Expiration Date:** 31 OCT 2026**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
Anti-rat IL-13 Antibody	D0082147	2-8°C	28 FEB 2027
Anti-rat TNF- α Antibody	D0082148	2-8°C	28 FEB 2027
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080437	2-8°C	30 JUN 2030
Proinflammatory Panel 1 (rat) Plate	Z0048027	2-8°C	28 FEB 2027
Rat IFN- γ Antibody	D0082353	2-8°C	29 FEB 2028
Rat IL-10 Antibody	D0082146	2-8°C	30 JUN 2027
Rat IL-1 β Antibody	D0082140	2-8°C	30 JUN 2027
Rat IL-2 Antibody	D0082141	2-8°C	30 JUN 2027
Rat IL-4 Antibody	D0082142	2-8°C	30 JUN 2027
Rat IL-5 Antibody	D0082143	2-8°C	30 JUN 2027
Rat IL-6 Antibody	D0082144	2-8°C	30 JUN 2027
Rat KC/GRO Antibody	D0082145	2-8°C	31 JUL 2027
Diluent 3	M0030140	$\leq -10^{\circ}\text{C}$	30 SEP 2027
Diluent 65	M0650002	$\leq -10^{\circ}\text{C}$	31 OCT 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 (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%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
2	IL-1 β	80% - 120%	< 20%	70% - 130%	< 25%	70% - 130%	< 25%	70% - 130%	< 25%
3	IL-2	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
4	IL-4	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
5	IL-5	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
6	IL-6	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
7	KC/GRO	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
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

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	1.21	75% - 125%	< 25%	292	75% - 125%	< 25%
2	IL-1 β	2.66	75% - 125%	< 25%	1230	75% - 125%	< 25%
3	IL-2	1.07	75% - 125%	< 25%	455	75% - 125%	< 25%
4	IL-4	3.62	75% - 125%	< 25%	357	75% - 125%	< 25%
5	IL-5	1.86	75% - 125%	< 25%	305	75% - 125%	< 25%
6	IL-6	188	75% - 125%	< 25%	54600	75% - 125%	< 25%
7	KC/GRO	2.23	75% - 125%	< 25%	604	75% - 125%	< 25%
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%



CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	453	pg/mL
2	IL-1 β	1690	pg/mL
3	IL-2	768	pg/mL
4	IL-4	573	pg/mL
5	IL-5	476	pg/mL
6	IL-6	78900	pg/mL
7	KC/GRO	1060	pg/mL
8	IL-10	7590	pg/mL
9	IL-13	1500	pg/mL
10	TNF- α	1260	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		10 SEP 2025