



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 2 (rat) Kit
Lot Number: K0082551
Expiration Date: 30 SEP 2025
Catalog Number: K15059-Series; K153QP-Series; K153QU-Series; K153QO-Series; K153QX-Series; K153QR-Series; K153QS-Series; K153QT-Series; K153OD-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat IFN- γ Antibody	D0082239	2-8°C	30 SEP 2027
Anti-rat IL-10 Antibody	D0082245	2-8°C	30 SEP 2027
Anti-rat IL-13 Antibody	D0082246	2-8°C	30 SEP 2027
Anti-rat IL-1 β Antibody	D0082241	2-8°C	30 SEP 2027
Anti-rat IL-4 Antibody	D0082150	2-8°C	31 MAR 2027
Anti-rat IL-5 Antibody	D0082242	2-8°C	30 SEP 2027
Anti-rat IL-6 Antibody	D0082243	2-8°C	30 SEP 2027
Anti-rat KC/GRO Antibody	D0082244	2-8°C	30 SEP 2027
Anti-rat TNF- α Antibody	D0082247	2-8°C	30 SEP 2027
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080386	2-8°C	28 FEB 2029
Proinflammatory Panel 2 (rat) Plate	Z0047974	2-8°C	30 SEP 2026
Diluent 40	M0400020	$\leq -10^{\circ}\text{C}$	28 FEB 2026
Diluent 42	M0420019	$\leq -10^{\circ}\text{C}$	30 SEP 2025
Blocker H	R0100024	Room Temperature	28 FEB 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%	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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	39.7	75% - 125%	< 25%	3750	75% - 125%	< 25%
2	IL-1 β	102	75% - 125%	< 25%	8100	75% - 125%	< 25%
4	IL-4	8	75% - 125%	< 25%	723	75% - 125%	< 25%
5	IL-5	82	75% - 125%	< 25%	3000	75% - 125%	< 25%
6	IL-6	96.9	75% - 125%	< 25%	8550	75% - 125%	< 25%
7	KC/GRO	21.7	75% - 125%	< 25%	728	75% - 125%	< 25%
8	IL-10	163	75% - 125%	< 25%	15700	75% - 125%	< 25%
9	IL-13	12.5	75% - 125%	< 25%	1080	75% - 125%	< 25%
10	TNF- α	9.1	75% - 125%	< 25%	793	75% - 125%	< 25%

Non-Specific Binding Specifications:

Spot	Name	Spec
1	IFN- γ	< 1.0%
2	IL-1 β	< 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	5230	pg/mL
2	IL-1 β	12700	pg/mL
4	IL-4	1110	pg/mL
5	IL-5	8440	pg/mL
6	IL-6	13900	pg/mL
7	KC/GRO	2700	pg/mL
8	IL-10	30000	pg/mL
9	IL-13	1560	pg/mL
10	TNF- α	1200	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		18 SEP 2024