

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 2 (rat) Kit
Kit Lot Number: K0082159
Expiration Date: 31 Jul 2024
Kit Catalog Numbers: K15059-Series; K153OD-Series; K153QO-Series; K153QP-Series; K153QR-Series; K153QS-Series; K153QT-Series; K153QU-Series; K153QX-Series

Precision and Accuracy 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%

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	8.00	75% - 125%	< 25%	723	75% - 125%	< 25%
5	IL-5	82.0	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	9.10	75% - 125%	< 25%	793	75% - 125%	< 25%

Non-Specific Binding Specifications:

Spot	Name	NSB 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%

CERTIFICATE OF ANALYSIS

Page 2 of 2

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IFN- γ	5240	pg/mL
2	IL-1 β	20500	pg/mL
4	IL-4	988	pg/mL
5	IL-5	9030	pg/mL
6	IL-6	13500	pg/mL
7	KC/GRO	2450	pg/mL
8	IL-10	27700	pg/mL
9	IL-13	1500	pg/mL
10	TNF- α	1150	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 2 (rat) Plate	Z0047782	2-8°C	30 Apr 2025
Anti-rat IFN- γ Antibody	D0081939	2-8°C	31 Mar 2026
Anti-rat IL-1 β Antibody	D0081941	2-8°C	31 Mar 2026
Anti-rat IL-4 Antibody	D0081940	2-8°C	31 Mar 2026
Anti-rat IL-5 Antibody	D0081942	2-8°C	31 Mar 2026
Anti-rat IL-6 Antibody	D0081943	2-8°C	31 Mar 2026
Anti-rat KC/GRO Antibody	D0081944	2-8°C	31 Mar 2026
Anti-rat IL-10 Antibody	D0081945	2-8°C	31 Mar 2026
Anti-rat IL-13 Antibody	D0081946	2-8°C	31 Mar 2026
Anti-rat TNF- α Antibody	D0081947	2-8°C	31 Mar 2026
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080337	2-8°C	30 Sep 2027
Blocker H	R0100022	Room Temperature	31 Oct 2024
Diluent 40	M0400018	$\leq -10^{\circ}\text{C}$	31 Jul 2024
Diluent 42	M0420016	$\leq -10^{\circ}\text{C}$	31 Jul 2024

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 DEC 2022