

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 2 (rat) Kit
Kit Lot Number: K0081905
Expiration Date: 31 Jul 2023
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

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IFN- γ	6030	pg/mL
2	IL-1 β	14700	pg/mL
4	IL-4	1110	pg/mL
5	IL-5	10400	pg/mL
6	IL-6	15100	pg/mL
7	KC/GRO	2610	pg/mL
8	IL-10	27000	pg/mL
9	IL-13	1170	pg/mL
10	TNF- α	1170	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 2 (rat) Plate	Z0047495	2-8°C	30 Apr 2024
Anti-rat IFN- γ Antibody	D0081793	2-8°C	31 Oct 2025
Anti-rat IL-1 β Antibody	D0081736	2-8°C	31 Jan 2025
Anti-rat IL-4 Antibody	D0081735	2-8°C	31 Jan 2025
Anti-rat IL-5 Antibody	D0081737	2-8°C	31 Mar 2025
Anti-rat IL-6 Antibody	D0081738	2-8°C	31 Jan 2025
Anti-rat KC/GRO Antibody	D0081739	2-8°C	28 Feb 2025
Anti-rat IL-10 Antibody	D0081740	2-8°C	31 Mar 2025
Anti-rat IL-13 Antibody	D0081741	2-8°C	28 Feb 2025
Anti-rat TNF- α Antibody	D0081742	2-8°C	31 Jan 2025
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080313	2-8°C	31 Mar 2027
Blocker H	R0100021	Room Temperature	30 Nov 2023
Diluent 42	M0420015	$\leq -10^\circ\text{C}$	31 Mar 2024
Diluent 40	M0400017	$\leq -10^\circ\text{C}$	31 Jul 2023

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		25 AUG 2022