



Extension Memo

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 2 (rat) Kit
Lot Number: K0082908
New Expiration Date: 31 AUG 2027
Catalog Number(s): K153QP-Series; K153QX-Series; K153QU-Series; K15059-Series; K153QO-Series;
 K153QR-Series; K153QS-Series; K153QT-Series; K153OD-Series

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 42	M0420021	31 MAR 2027	31 AUG 2027
Diluent 40	M0400022	30 NOV 2026	31 AUG 2027
Blocker H	R0100025	31 AUG 2026	31 AUG 2027

Function	Name	Signature	Date
Quality	Michelle Likiam		02 Jun 2026

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 2 (rat) Kit
Lot Number: K0082908
Expiration Date: 31 AUG 2026
Catalog Number: K153QP-Series; K153QX-Series; K153QU-Series; K15059-Series; K153QO-Series;
 K153QR-Series; K153QS-Series; K153QT-Series; K153OD-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat IFN- γ Antibody	D0082385	2-8°C	31 JUL 2028
Anti-rat IL-10 Antibody	D0082391	2-8°C	31 JUL 2028
Anti-rat IL-13 Antibody	D0082392	2-8°C	31 MAY 2028
Anti-rat IL-1 β Antibody	D0082387	2-8°C	31 MAY 2028
Anti-rat IL-4 Antibody	D0082386	2-8°C	31 MAY 2028
Anti-rat IL-5 Antibody	D0082388	2-8°C	30 JUN 2028
Anti-rat IL-6 Antibody	D0082389	2-8°C	31 MAY 2028
Anti-rat KC/GRO Antibody	D0082390	2-8°C	30 JUN 2028
Anti-rat TNF- α Antibody	D0082393	2-8°C	30 NOV 2028
Proinflammatory Panel 1(rat) Calibrator Blend	A0080413	2-8°C	30 APR 2030
Proinflammatory Panel 2 (rat) 96-Well 10-Spot SECTOR Plate	Z0048241	2-8°C	31 AUG 2028
Diluent 40	M0400022	$\leq -10^{\circ}\text{C}$	30 NOV 2026
Diluent 42	M0420021	$\leq -10^{\circ}\text{C}$	31 MAR 2027
Blocker H	R0100025	Room Temperature	31 AUG 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	6300	pg/mL
2	IL-1 β	11900	pg/mL
4	IL-4	1050	pg/mL
5	IL-5	9540	pg/mL
6	IL-6	13400	pg/mL
7	KC/GRO	2950	pg/mL
8	IL-10	23900	pg/mL
9	IL-13	1670	pg/mL
10	TNF- α	1330	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		02 JUN 2026