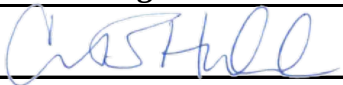




Extension Memo

Product Description: Diluent 65
Lot Number: M0650002
New Expiration Date: 31 OCT 2027
Catalog Number(s): R50CJ-Series

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

Function	Name	Signature	Date
Quality	Carissa B. Hubbard		10 Jul 2026

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Proinflammatory Panel 1 (rat) Kit, SECTOR**Lot Number:** K0082969**Expiration Date:** 28 FEB 2027**Catalog Number:** K15294-Series, K153J2-Series, K153K2-Series, K153L2-Series, K153M2-Series, K153N2-Series, K153P2-Series, K153Q2-Series, K153R2-Series, K153S2-Series, K153B2-Series, K15325-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 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) 96-Well 10-Spot SECTOR Plate	Z0048027	2-8°C	28 FEB 2027
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080380	2-8°C	31 MAR 2029
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 2027

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%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	456	pg/mL
2	IL-1 β	2180	pg/mL
3	IL-2	758	pg/mL
4	IL-4	544	pg/mL
5	IL-5	478	pg/mL
6	IL-6	84500	pg/mL
7	KC/GRO	1110	pg/mL
8	IL-10	6610	pg/mL
9	IL-13	1540	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	Carissa B. Hubbard	Quality		16 JUL 2026