



RFE MEMO

Page 1 of 1

Product Description(s): Diluent 42

Catalog Number(s) / Size(s): R50AK-Series

Lot Number: M0420010

Based on additional testing, the original date of expiration for this diluent has been extended by 12 months. The expiration date of M0420010 has been extended to 31 JUL 2021.

Function	Name	Signature	Date
Quality	Karen Hamilla		07 APR 2020

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 2 (Rat) Kit
Kit Lot Number: K0081372
Expiration Date: 31 Jul 2021
Kit Catalog Numbers: K15059-Series, K153OD-Series, K153QO-Series, K153QR-Series, K153QS-Series, K153QT-Series

Precision and Accuracy Specifications:

		Calibrator		Proinflammatory Panel 1 (rat) Control 1		Proinflammatory Panel 1 (rat) Control 2		Proinflammatory Panel 1 (rat) Control 3	
Spot	Name	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%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	102	75% - 125%	< 25%	8100	75% - 125%	< 25%
2	IL-1 β	8.00	75% - 125%	< 25%	723	75% - 125%	< 25%
4	IL-4	82.0	75% - 125%	< 25%	3000	75% - 125%	< 25%
5	IL-5	96.9	75% - 125%	< 25%	8550	75% - 125%	< 25%
6	IL-6	21.7	75% - 125%	< 25%	728	75% - 125%	< 25%
7	KC/GRO	163	75% - 125%	< 25%	15700	75% - 125%	< 25%
8	IL-10	12.5	75% - 125%	< 25%	1080	75% - 125%	< 25%
9	IL-13	9.10	75% - 125%	< 25%	793	75% - 125%	< 25%
10	TNF- α	0	0	0	0	0	0

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	6100	pg/mL
2	IL-1 β	19500	pg/mL
4	IL-4	1100	pg/mL
5	IL-5	9440	pg/mL
6	IL-6	9830	pg/mL
7	KC/GRO	2680	pg/mL
8	IL-10	31400	pg/mL
9	IL-13	1550	pg/mL
10	TNF- α	1150	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 2 (Rat) Plate	Z0047221	2-8°C	30 Apr 2022
Anti-rat IFN- γ Antibody	D0081405	2-8°C	30 Jun 2023
Anti-rat IL-1 β Antibody	D0081407	2-8°C	31 May 2023
Anti-rat IL-4 Antibody	D0081406	2-8°C	30 Apr 2023
Anti-rat IL-5 Antibody	D0081408	2-8°C	30 Jun 2023
Anti-rat IL-6 Antibody	D0081350	2-8°C	31 May 2023
Anti-rat KC/GRO Antibody	D0081410	2-8°C	30 Apr 2023
Anti-rat IL-10 Antibody	D0081411	2-8°C	30 Apr 2023
Anti-rat IL-13 Antibody	D0081412	2-8°C	30 Apr 2023
Anti-rat TNF- α Antibody	D0081413	2-8°C	30 Apr 2023
Proinflammatory Panel 1 (Rat) Calibrator Blend	A0080248	2-8°C	31 Jan 2025
Blocker H	R0100018	Room Temperature	31 Mar 2022
Diluent 42	M0420010	$\leq -10^{\circ}\text{C}$	31 Jul 2021
Diluent 40	M0400014	$\leq -10^{\circ}\text{C}$	31 Aug 2021

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		01 FEB 2021