



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Cytokine Panel 1 (human) Kit
Catalog Number(s) / Size(s): K15050-Series; K15057-Series; K151RC-Series; K151RG-Series;
K156RC-Series; K156RG-Series; K151RE-Series; K156RE-Series
Lot Number: K0081762

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081762 has been extended to 28 FEB 2023.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 43	M0430034	28 FEB 2022	28 FEB 2023

Function	Name	Signature	Date
Quality	Karen Hamilla		13 DEC 2021

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 Cytokine Panel 1 (human) Kit
Kit Lot Number: K0081762
Expiration Date: 28 Feb 2022
Kit Catalog Numbers: K15050-Series, K15057-Series, K151RC-Series, K151RG-Series, K156RC-Series, K156RG-Series, K151RE-Series, K156RE-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	GM-CSF	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	IL-1 α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	IL-5	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	IL-7	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	IL-12/IL-23p40	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	IL-15	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	IL-16	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	IL-17A	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	TNF- β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	VEGF	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	GM-CSF	0.842	80% - 120%	< 20%	750	80% - 120%	< 20%
2	IL-1 α	2.85	80% - 120%	< 20%	278	80% - 120%	< 20%
3	IL-5	4.41	80% - 120%	< 20%	562	80% - 120%	< 20%
4	IL-7	0.851	80% - 120%	< 20%	563	80% - 120%	< 20%
5	IL-12/IL-23p40	1.32	80% - 120%	< 20%	2250	80% - 120%	< 20%
6	IL-15	0.774	80% - 120%	< 20%	525	80% - 120%	< 20%
7	IL-16	19.1	80% - 120%	< 20%	1870	80% - 120%	< 20%
8	IL-17A	3.19	80% - 120%	< 20%	3650	80% - 120%	< 20%
9	TNF- β	0.465	80% - 120%	< 20%	458	80% - 120%	< 20%
10	VEGF	7.70	80% - 120%	< 20%	562	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	NSB Spec
1	GM-CSF	< 1.0%
2	IL-1 α	< 1.0%
3	IL-5	< 1.0%
4	IL-7	< 1.0%
5	IL-12/IL-23p40	< 1.0%
6	IL-15	< 1.0%
7	IL-16	< 1.0%
8	IL-17A	< 1.0%
9	TNF- β	< 1.0%
10	VEGF	< 1.0%

CERTIFICATE OF ANALYSIS

Page 2 of 2

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	GM-CSF	1200	pg/mL
2	IL-1 α	438	pg/mL
3	IL-5	853	pg/mL
4	IL-7	766	pg/mL
5	IL-12/IL-23p40	3270	pg/mL
6	IL-15	714	pg/mL
7	IL-16	2930	pg/mL
8	IL-17A	6710	pg/mL
9	TNF- β	792	pg/mL
10	VEGF	1150	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Plate	Z0047512	2-8°C	30 Nov 2023
Anti-hu GM-CSF Antibody	D0081654	2-8°C	31 Aug 2024
Anti-hu IL-1 α Antibody	D0081644	2-8°C	30 Sep 2024
Anti-hu IL-5 Antibody	D0081645	2-8°C	30 Sep 2024
Anti-hu IL-7 Antibody	D0081530	2-8°C	30 Jun 2024
Anti-hu IL-12/23p40 Antibody	D0081671	2-8°C	30 Nov 2024
Anti-hu IL-15 Antibody	D0081655	2-8°C	30 Sep 2024
Anti-hu IL-16 Antibody	D0081611	2-8°C	30 Jun 2024
Anti-hu IL-17A Antibody	D0081656	2-8°C	31 Aug 2024
Anti-hu TNF- β Antibody	D0081647	2-8°C	30 Nov 2024
Anti-hu VEGF Antibody	D0081574	2-8°C	31 Mar 2024
Cytokine Panel 1 (human) Calibrator Blend	A0080283	2-8°C	31 Mar 2026
Diluent 43	M0430034	$\leq -10^\circ\text{C}$	28 Feb 2022
Diluent 3	M0030104	$\leq -10^\circ\text{C}$	30 Jun 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		02 AUG 2021