



EXTENSION MEMO

Page 1 of 1

Product Description: V-PLEX® and V-PLEX plus Human GM-CSF Kit


Catalog Number: K156RIG-Series; K151RID-Series

Lot Number: K0082357

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082357 has been extended to 30 NOV 2024.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human GM-CSF Plate	Z0047553	30 Nov 2023	30 NOV 2024
Anti-hu GM-CSF Antibody	D0081654	31 Aug 2024	30 NOV 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		14 SEP 2023

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

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human GM-CSF Kit

Kit Lot Number: K0082357

Expiration Date: 30 Nov 2023

Kit Catalog Numbers: K151RI-Series; K156RI-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human GM-CSF Plate	Z0047553	2-8°C	30 Nov 2023
Anti-hu GM-CSF Antibody	D0081654	2-8°C	31 Aug 2024
Cytokine Panel 1(human) Calibrator Blend	A0080300	2-8°C	31 Jul 2026
Diluent 3	M0030120	≤ -10°C	31 Jul 2025
Diluent 43	M0430050	≤ -10°C	31 Mar 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%


Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	GM-CSF	0.842	80% - 120%	< 20%	750	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	GM-CSF	1290	pg/mL

Dilution to Top of Curve = 1X

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 Sep 2023