

CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Kit Lot Number: K0082358

Expiration Date: 30 Apr 2024

Kit Catalog Numbers: K151RI-Series; K156RI-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human GM-CSF Plate	Z0047623	2-8°C	30 Apr 2024
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