



CERTIFICATE OF ANALYSIS

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

Lot Number: K0082673

Expiration Date: 30 SEP 2026

Catalog Number: K151RI-Series; K156RI-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu GM-CSF Antibody	D0082095	2-8°C	31 DEC 2026
Cytokine Panel 1 (human) Calibrator Blend	A0080383	2-8°C	31 DEC 2028
Human GM-CSF 96-Well Small Spot SECTOR Plate	Z0048098	2-8°C	31 AUG 2027
Diluent 3	M0030136	≤ -10°C	31 MAY 2027
Diluent 43	M0430056	≤ -10°C	30 SEP 2026

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

Precision, Accuracy, and Sensitivity 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	1110	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	Karen Hamilla	Quality		07 APR 2025