



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX® Human GM-CSF Kit

Catalog Number: K151F3S-Series

Lot Number: K00S0160

Expiration Date: 31 JAN 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human GM-CSF Antibody	B00T0169	2-8°C	31 JAN 2027
TURBO-BOOST™ Human GM-CSF Antibody	D00T0169	2-8°C	28 FEB 2027
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
Blocker S1	R0010832	≤ -10°C	31 OCT 2025
S-PLEX Enhance E1	Y0550012	≤ -10°C	31 AUG 2025
S-PLEX Enhance E2	Y0560010	≤ -10°C	31 JAN 2025
S-PLEX Enhance E3	Y0570014	≤ -70°C	31 OCT 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580018	≤ -70°C	31 OCT 2025
S-PLEX Detect D2	Y0590014	≤ -70°C	31 OCT 2025
Human GM-CSF Calibrator	A00S0043	2-8°C	31 DEC 2025
Diluent 43	M0430053	≤ -10°C	31 JAN 2026
Diluent 3	M0030123	≤ -10°C	31 JAN 2026

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

Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator	
	Recovery	CV
GM-CSF	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
GM-CSF	0.550 - 8.80	8.80	fg/mL	80% - 120%	< 20%	2.01	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
GM-CSF	50X Calibrator	1350000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		22 NOV 2023