



## CERTIFICATE OF ANALYSIS

**Product Description:** S-PLEX<sup>®</sup> Human GM-CSF Kit  
**Lot Number:** K00S0242  
**Expiration Date:** 31 AUG 2025  
**Catalog Number:** K151F3S-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human GM-CSF Antibody	B00T0169	2-8°C	31 JAN 2027
Human GM-CSF Calibrator	A00S0043	2-8°C	31 DEC 2025
TURBO-BOOST Human GM-CSF Antibody	D00T0169	2-8°C	28 FEB 2027
Blocker S1	R0010832	≤ -10°C	31 OCT 2025
Diluent 3	M0030123	≤ -10°C	31 JAN 2026
Diluent 43	M0430053	≤ -10°C	31 JAN 2026
S-PLEX Enhance E1	Y0550012	≤ -10°C	31 AUG 2025
S-PLEX Enhance E2	Y0560019	≤ -10°C	30 NOV 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> Reagent	Y0580018	≤ -70°C	31 OCT 2025
S-PLEX Detect D2	Y0590014	≤ -70°C	31 OCT 2025
S-PLEX Enhance E3	Y0570014	≤ -70°C	31 OCT 2025

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

**Precision, Accuracy, and Sensitivity Specifications:**

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

		LLOQ				LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
GM-CSF	0.550 - 8.80	8.80	fg/mL	80% - 120%	< 20%	2.34	fg/mL

**Stock Concentration of Calibrator:**

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

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		22 NOV 2024