

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human GM-CSF Kit
Lot Number: K00S0237
Expiration Date: 31 DEC 2025
Catalog Number: K151F3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human GM-CSF Antibody	B00T0237	2-8°C	30 JUN 2028
Human GM-CSF Calibrator	A00S0043	2-8°C	31 DEC 2025
TURBO-BOOST Human GM-CSF Antibody	D00T0237	2-8°C	31 MAY 2028
Blocker S1	R0140011	≤ -10°C	28 FEB 2027
Diluent 3	M0030130	≤ -10°C	31 JAN 2027
Diluent 43	M0430056	≤ -10°C	30 SEP 2026
S-PLEX Enhance E1	Y0550020	≤ -10°C	30 APR 2027
S-PLEX Enhance E2	Y0560020	≤ -10°C	31 MAR 2027
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580029	≤ -70°C	30 APR 2027
S-PLEX Detect D2	Y0590022	≤ -70°C	30 APR 2027
S-PLEX Enhance E3	Y0570023	≤ -70°C	30 APR 2027

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

Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
GM-CSF	0.550 - 8.80	8.80	fg/mL	80% - 120%	< 20%	2.81	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
Review/Approval	Karen Hamilla	Quality		20 FEB 2025