



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151F3S-Series

Lot Number: K00S0091

Expiration Date: 30 APR 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human GM-CSF Antibody	D00T0091	2-8°C	31 MAR 2026
Biotin Human GM-CSF Antibody	B00T0091	2-8°C	28 FEB 2026
Human GM-CSF Calibrator	A00S0043	2-8°C	31 DEC 2025
Diluent 43	M0430047	≤ -10°C	31 MAY 2024
Diluent 3	M0030116	≤ -10°C	30 NOV 2024
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0010811	≤ -10°C	28 FEB 2025
S-PLEX Enhance E1	Y0550009	≤ -10°C	28 FEB 2025
S-PLEX Enhance E2	Y0560011	≤ -10°C	31 MAR 2025
S-PLEX Enhance E3	Y0570011	≤ -70°C	31 MAR 2025
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580015	≤ -70°C	30 APR 2025
S-PLEX Detect D2	Y0590011	≤ -70°C	31 MAR 2025

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	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
GM-CSF	0.550 – 8.80	8.80	fg/mL	80% - 120%	< 20%	1.66	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		30 DEC 2022