

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human G-CSF Kit
Catalog Number: K151K3S-Series
Lot Number: K00S0187
Expiration Date: 31 JUL 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human G-CSF Antibody	B00T0121	2-8°C	31 MAY 2026
Human G-CSF Calibrator	A00S0015	2-8°C	31 JUL 2024
TURBO-BOOST™ Human G-CSF Antibody	D00T0121	2-8°C	31 MAY 2026
Blocker S1	R0010832	≤ -10°C	31 OCT 2025
Diluent 3	M0030125	≤ -10°C	31 MAY 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026
S-PLEX Enhance E1	Y0550013	≤ -10°C	30 NOV 2025
S-PLEX Enhance E2	Y0560015	≤ -10°C	28 FEB 2026
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580019	≤ -70°C	31 DEC 2025
S-PLEX Detect D2	Y0590015	≤ -70°C	31 JAN 2026
S-PLEX Enhance E3	Y0570015	≤ -70°C	30 NOV 2025

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	G-CSF	75% - 125%	< 25%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
G-CSF	27.5 - 440	55.0	fg/mL	75% - 125%	< 25%	42.5	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
G-CSF	50X Calibrator	27000000	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		28 FEB 2024