

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human G-CSF Kit
Lot Number: K00S0190
Expiration Date: 31 OCT 2025
Catalog Number: K151K3S-Series; K156K3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human G-CSF Antibody	B00T0199	2-8°C	31 OCT 2027
TURBO-BOOST™ Human G-CSF Antibody	D00T0199	2-8°C	31 OCT 2027
Human G-CSF Calibrator	A00S0105	2-8°C	31 MAR 2029
Diluent 3	M0030128	≤ -10°C	31 OCT 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026
Blocker S1	R0010832	≤ -10°C	31 OCT 2025
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580023	≤ -70°C	31 JUL 2026
S-PLEX Detect D2	Y0590017	≤ -70°C	30 JUN 2026
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560018	≤ -10°C	31 AUG 2026
S-PLEX Enhance E3	Y0570018	≤ -70°C	30 JUN 2026

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	110	fg/mL	75% - 125%	< 25%	66.6	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
G-CSF	50X Calibrator	26450000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		04 SEP 2024