



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human G-CSF Kit  
**Catalog Number:** K151K3S-Series; K156K3S-Series  
**Lot Number:** K00S0113  
**Expiration Date:** 30 APR 2024

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human G-CSF Antibody	D00T0072	2-8°C	30 NOV 2025
Biotin Human G-CSF Antibody	B00T0072	2-8°C	30 NOV 2025
Human G-CSF Calibrator	A00S0015	2-8°C	31 JUL 2024
Diluent 3	M0030116	≤ -10°C	30 NOV 2024
Diluent 43	M0430047	≤ -10°C	31 MAY 2024
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0010768	≤ -10°C	31 MAY 2024
S-PLEX Enhance E1	Y0550008	≤ -10°C	30 NOV 2024
S-PLEX Enhance E2	Y0560010	≤ -10°C	31 JAN 2025
S-PLEX Enhance E3	Y0570010	≤ -70°C	31 JAN 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580014	≤ -70°C	28 FEB 2025
S-PLEX Detect D2	Y0590010	≤ -70°C	28 FEB 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
G-CSF	75% - 125%	< 25%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
G-CSF	27.5 - 440	110	fg/mL	75% - 125%	< 25%	44.0	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
G-CSF	50X Calibrator	27000000	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		10 NOV 2022