



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human G-CSF Kit
Catalog Number: K151K3S-Series; K156K3S-Series
Lot Number: K00S0120
Expiration Date: 29 FEB 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human G-CSF Antibody	D00T0121	2-8°C	31 MAY 2026
Biotin Human G-CSF Antibody	B00T0121	2-8°C	31 MAY 2026
Human G-CSF Calibrator	A00S0015	2-8°C	31 JUL 2024
Diluent 3	M0030111	≤ -10°C	31 MAY 2024
Diluent 43	M0430046	≤ -10°C	29 FEB 2024
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0010811	≤ -10°C	28 FEB 2025
S-PLEX Enhance E1	Y0550010	≤ -10°C	31 MAY 2025
S-PLEX Enhance E2	Y0560008	≤ -10°C	30 APR 2024
S-PLEX Enhance E3	Y0570011	≤ -70°C	30 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
G-CSF	75% - 125%	< 25%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
G-CSF	27.5 - 440	55.0	fg/mL	75% - 125%	< 25%	41.0	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		13 JAN 2023