



### 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:** K00S0021  
**Expiration Date:** 31 DEC 2021

#### Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
G-CSF	75% - 125%	< 25%

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

LLOD		
Name	Observed	Units
G-CSF	58.8	fg/mL

#### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
G-CSF	50X Calibrator	27000000	fg/mL

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human G-CSF Antibody	D00T0012A	2-8°C	31 Oct 2022
Biotin Human G-CSF Antibody	B00T0012A	2-8°C	31 Oct 2022
Human G-CSF Calibrator	A00S0015	2-8°C	31 Jul 2024
Diluent 43	M0430033	≤ -10°C	31 Jan 2022
Diluent 3	M0030091	≤ -10°C	31 Jan 2022
S-PLEX Coating Reagent C1	288327	≤ -70°C	31 May 2022
Blocker S1	R0140002	≤ -10°C	31 Dec 2021
S-PLEX Enhance E1	Y0550002	≤ -10°C	31 Dec 2021
S-PLEX Enhance E2	Y0560002	≤ -10°C	31 Dec 2021
S-PLEX Enhance E3	Y0570002	≤ -70°C	31 Dec 2021
S-PLEX Detect D1	Y0580002	≤ -70°C	31 Jan 2022
S-PLEX Detect D2	Y0590003	≤ -70°C	31 Aug 2022

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		22 JUN 2020