



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human G-CSF Kit

Catalog Number: K151K3S-Series; K156K3S-Series

Lot Number: K00S0013

Expiration Date: 31 OCT 2020

Precision, Accuracy, and Sensitivity Specifications:

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

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

Name	LLOD	
	Observed	Units
G-CSF	44.1	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
S-PLEX [®] 96-Well SECTOR Plate	Z00S0003	2-8°C	31 DEC 2021
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	Y0590002	≤ -70°C	31 OCT 2020

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
Review/Approval	Karen Hamilla	Quality		29 JUL 2020