



Extension Memo

Page 1 of 1

Product Description: S-PLEX Human G-CSF Kit
Catalog Number(s): K151K3S-Series, K156K3S-Series
Lot Number: K00S0190
New Expiration Date: 31 Aug 2026

Based on additional testing, the original date of expiration for this kit has been extended by 10 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 43	M0430055	30 Apr 2026	31 Aug 2026
Blocker S1	R0010832	31 Oct 2025	31 Aug 2026
S-PLEX Coating Reagent C1	446866	31 Mar 2026	31 Aug 2026
S-PLEX Detect D1 with TURBO-TAG reagent	Y0580023	31 Jul 2026	31 Aug 2026
S-PLEX Enhance E3	Y0570018	30 Jun 2026	31 Aug 2026
S-PLEX Detect D2	Y0590017	30 Jun 2026	31 Aug 2026
S-PLEX Enhance E1	Y0550016	31 Jul 2026	31 Aug 2026

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		11 Aug 2025



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:

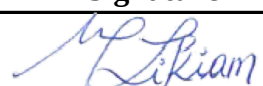
Spot Number	Name	Calibrator	
		Recovery	CV
1	G-CSF	75% - 125%	< 25%

Name	LLOQ					LLOD	
	Range Tested	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