

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human G-CSF Kit
Lot Number: K00S0292
Expiration Date: 30 SEP 2026
Catalog Number: K151K3S-Series; K156K3S-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human G-CSF Antibody	B00T0270	2-8°C	31 MAR 2029
Human G-CSF Calibrator	A00S0105	2-8°C	31 MAR 2029
TURBO-BOOST Human G-CSF Antibody	D00T0270	2-8°C	30 JUN 2029
Blocker S1	R0140017	≤ -10°C	29 FEB 2028
Diluent 3	M0030136	≤ -10°C	31 MAY 2027
Diluent 43	M0430056	≤ -10°C	30 SEP 2026
S-PLEX Enhance E1	Y0550027	≤ -10°C	29 FEB 2028
S-PLEX Enhance E2	Y0560026	≤ -10°C	29 FEB 2028
S-PLEX Coating Reagent C1	521329	≤ -70°C	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580035	≤ -70°C	31 JAN 2028
S-PLEX Detect D2	Y0590027	≤ -70°C	31 JAN 2028
S-PLEX Enhance E3	Y0570029	≤ -70°C	29 FEB 2028

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	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
G-CSF	27.5 - 440	55.0	fg/mL	75% - 125%	< 25%	35.0	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
G-CSF	50X Calibrator	26450000	fg/mL

This lot meets MSD's non-specific binding specifications.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 MAR 2026