

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX Human IFN- $\gamma$  Kit  
**Lot Number:** K00S0199  
**Expiration Date:** 31 JUL 2025  
**Catalog Number:** K151X9S-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- $\gamma$ Antibody	B00T0085	2-8°C	30 SEP 2025
Human IFN- $\gamma$ Calibrator	A00S0077	2-8°C	31 OCT 2027
TURBO-BOOST Human IFN- $\gamma$ Antibody	D00T0085	2-8°C	31 DEC 2025
Blocker S1	R0010831	$\leq -10^{\circ}\text{C}$	31 JUL 2025
Diluent 58	M0580011	$\leq -10^{\circ}\text{C}$	31 JUL 2026
S-PLEX Enhance E1	Y0550014	$\leq -10^{\circ}\text{C}$	28 FEB 2026
S-PLEX Enhance E2	Y0560015	$\leq -10^{\circ}\text{C}$	28 FEB 2026
S-PLEX Coating Reagent C1	417723	$\leq -70^{\circ}\text{C}$	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG <sup>TM</sup> Reagent	Y0580017	$\leq -70^{\circ}\text{C}$	31 AUG 2025
S-PLEX Detect D2	Y0590015	$\leq -70^{\circ}\text{C}$	31 JAN 2026
S-PLEX Enhance E3	Y0570013	$\leq -70^{\circ}\text{C}$	31 JUL 2025

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	IFN- $\gamma$	80% - 120%	< 20%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- $\gamma$	3.96 - 63.4	15.9	fg/mL	80% - 120%	< 20%	3.48	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
IFN- $\gamma$	1X Calibrator	32600	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		12 JUN 2024