

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human IFN- γ Kit
Catalog Number: K151X9S-Series
Lot Number: K00S0179
Expiration Date: 31 OCT 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- γ Antibody	B00T0184	2-8°C	30 JUN 2027
Human IFN- γ Calibrator	A00S0077	2-8°C	31 OCT 2027
TURBO-BOOST [™] Human IFN- γ Antibody	D00T0184	2-8°C	31 JUL 2027
Blocker S1	R0140007	$\leq -10^{\circ}\text{C}$	30 JUN 2026
Diluent 58	M0580011	$\leq -10^{\circ}\text{C}$	31 JUL 2026
S-PLEX Enhance E1	Y0550015	$\leq -10^{\circ}\text{C}$	31 MAY 2026
S-PLEX Enhance E2	Y0560015	$\leq -10^{\circ}\text{C}$	28 FEB 2026
S-PLEX Coating Reagent C1	446866	$\leq -70^{\circ}\text{C}$	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580018	$\leq -70^{\circ}\text{C}$	31 OCT 2025
S-PLEX Detect D2	Y0590017	$\leq -70^{\circ}\text{C}$	30 JUN 2026
S-PLEX Enhance E3	Y0570017	$\leq -70^{\circ}\text{C}$	31 MAR 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	IFN- γ	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IFN- γ	3.96 - 63.4	31.7	fg/mL	80% - 120%	< 20%	4.19	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- γ	1X Calibrator	32600	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 APR 2024