



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IFN- $\gamma$  Kit

**Catalog Number:** K151X9S - Series

**Lot Number:** K00S0105

**Expiration Date:** 31 MAR 2025

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human IFN- $\gamma$ Antibody	D00T0110	2-8°C	30 JUN 2026
Biotin Human IFN- $\gamma$ Antibody	B00T0110	2-8°C	31 MAR 2026
Human IFN- $\gamma$ Calibrator	A00S0077	2-8°C	31 OCT 2027
Diluent 58	M0580007	$\leq -10^{\circ}\text{C}$	30 APR 2025
S-PLEX Coating Reagent C1	417723	$\leq -70^{\circ}\text{C}$	31 AUG 2025
Blocker S1	R0010824	$\leq -10^{\circ}\text{C}$	30 JUN 2025
S-PLEX Enhance E1	Y0550011	$\leq -10^{\circ}\text{C}$	31 JUL 2025
S-PLEX Enhance E2	Y0560011	$\leq -10^{\circ}\text{C}$	31 MAR 2025
S-PLEX Enhance E3	Y0570012	$\leq -70^{\circ}\text{C}$	31 MAY 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580015	$\leq -70^{\circ}\text{C}$	30 APR 2025
S-PLEX Detect D2	Y0590011	$\leq -70^{\circ}\text{C}$	31 MAR 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

### Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator	
	Recovery	CV
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%	2.89	fg/mL

### Stock Concentration of Calibrator:

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 MAR 2023