



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IFN- γ Kit
Lot Number: K00S0219
Expiration Date: 31 JUL 2026
Catalog Number: K151X9S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- γ Antibody	B00T0226	2-8°C	31 DEC 2027
Human IFN- γ Calibrator	A00S0117	2-8°C	30 NOV 2029
TURBO-BOOST Human IFN- γ Antibody	D00T0226	2-8°C	29 FEB 2028
Blocker S1	R0140008	$\leq -10^{\circ}\text{C}$	31 JUL 2026
Diluent 58	M0580014	$\leq -10^{\circ}\text{C}$	31 MAY 2027
S-PLEX Enhance E1	Y0550020	$\leq -10^{\circ}\text{C}$	30 APR 2027
S-PLEX Enhance E2	Y0560020	$\leq -10^{\circ}\text{C}$	31 MAR 2027
S-PLEX Coating Reagent C1	516221	$\leq -70^{\circ}\text{C}$	31 OCT 2027
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580030	$\leq -70^{\circ}\text{C}$	30 APR 2027
S-PLEX Detect D2	Y0590023	$\leq -70^{\circ}\text{C}$	31 JUL 2027
S-PLEX Enhance E3	Y0570023	$\leq -70^{\circ}\text{C}$	30 APR 2027

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	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- γ	3.96 - 63.4	7.93	fg/mL	80% - 120%	< 20%	3.91	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- γ	1X Calibrator	33400	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		21 FEB 2025