

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human IFN- β Kit
Lot Number: K00S0212
Expiration Date: 31 JUL 2025
Catalog Number: K151ADRS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- β Antibody	B00T0137	2-8°C	31 AUG 2026
Human IFN- β Calibrator	A00S0101	2-8°C	31 DEC 2028
TURBO-BOOST Human IFN- β Antibody	D00T0137	2-8°C	31 AUG 2026
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	Y0550011	$\leq -10^{\circ}\text{C}$	31 JUL 2025
S-PLEX Enhance E2	Y0560013	$\leq -10^{\circ}\text{C}$	31 AUG 2025
S-PLEX Coating Reagent C1	417723	$\leq -70^{\circ}\text{C}$	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG TM reagent	Y0580017	$\leq -70^{\circ}\text{C}$	31 AUG 2025
S-PLEX Detect D2	Y0590013	$\leq -70^{\circ}\text{C}$	31 JUL 2025
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- β	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IFN- β	16.0 - 255	31.9	fg/mL	80% - 120%	< 20%	7.21	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- β	1X Calibrator	153000	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		03 JUN 2024