

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human IFN- β Kit
Lot Number: K00S0246
Expiration Date: 31 OCT 2026
Catalog Number: K151ADRS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- β Antibody	B00T0200	2-8°C	31 OCT 2027
Human IFN- β Calibrator	A00S0108	2-8°C	31 MAY 2029
TURBO-BOOST Human IFN- β Antibody	D00T0200	2-8°C	31 AUG 2027
Blocker S1	R0140009	$\leq -10^{\circ}\text{C}$	31 OCT 2026
Diluent 58	M0580012	$\leq -10^{\circ}\text{C}$	31 OCT 2026
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	476750	$\leq -70^{\circ}\text{C}$	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG [™] Reagent	Y0580029	$\leq -70^{\circ}\text{C}$	30 APR 2027
S-PLEX Detect D2	Y0590022	$\leq -70^{\circ}\text{C}$	30 APR 2027
S-PLEX Enhance E3	Y0570022	$\leq -70^{\circ}\text{C}$	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- β	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- β	16.0 - 255	31.9	fg/mL	80% - 120%	< 20%	10.1	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- β	1X Calibrator	155000	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 JAN 2025