



## CERTIFICATE OF ANALYSIS

**Product Description:** S-PLEX Human IFN- $\beta$  Kit  
**Catalog Number:** K151ADRS-Series  
**Lot Number:** K00S0161  
**Expiration Date** 31 JUL 2025

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- $\beta$ Antibody	B00T0170	2-8°C	31 JAN 2027
Human IFN- $\beta$ Calibrator	A00S0081	2-8°C	29 FEB 2028
TURBO-BOOST™ Human IFN- $\beta$ Antibody	D00T0170	2-8°C	28 FEB 2027
Blocker S1	R0010831	$\leq -10^{\circ}\text{C}$	31 JUL 2025
Diluent 58	M0580009	$\leq -10^{\circ}\text{C}$	30 NOV 2025
S-PLEX Enhance E1	Y0550012	$\leq -10^{\circ}\text{C}$	31 AUG 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™ reagent	Y0580018	$\leq -70^{\circ}\text{C}$	31 OCT 2025
S-PLEX Detect D2	Y0590014	$\leq -70^{\circ}\text{C}$	31 OCT 2025
S-PLEX Enhance E3	Y0570014	$\leq -70^{\circ}\text{C}$	31 OCT 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- $\beta$	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- $\beta$	15.9 - 255	15.9	fg/mL	80% - 120%	< 20%	5.57	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
IFN- $\beta$	1X Calibrator	157000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		11 DEC 2023