

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human IFN- β Kit
Catalog Number: K151ADRS-Series
Lot Number: K00S0186
Expiration Date: 31 JUL 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- β Antibody	B00T0170	2-8°C	31 JAN 2027
Human IFN- β Calibrator	A00S0101	2-8°C	31 DEC 2028
TURBO-BOOST™ Human IFN- β Antibody	D00T0170	2-8°C	28 FEB 2027
Blocker S1	R0010831	≤ -10°C	31 JUL 2025
Diluent 58	M0580009	≤ -10°C	30 NOV 2025
S-PLEX Enhance E1	Y0550012	≤ -10°C	31 AUG 2025
S-PLEX Enhance E2	Y0560013	≤ -10°C	31 AUG 2025
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580018	≤ -70°C	31 OCT 2025
S-PLEX Detect D2	Y0590014	≤ -70°C	31 OCT 2025
S-PLEX Enhance E3	Y0570014	≤ -70°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- β	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- β	15.9 - 255	15.9	fg/mL	80% - 120%	< 20%	6.89	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		23 FEB 2024