



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IFN- β Kit
Lot Number: K00S0260
Expiration Date: 31 MAY 2027
Catalog Number: K151ADRS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- β Antibody	B00T0245	2-8°C	31 OCT 2028
TURBO-BOOST Human IFN- β Antibody	D00T0245	2-8°C	30 NOV 2028
Human IFN- β Calibrator	A00S0108	2-8°C	31 MAY 2029
Blocker S1	R0140016	$\leq -10^{\circ}\text{C}$	31 DEC 2027
Diluent 58	M0580014	$\leq -10^{\circ}\text{C}$	31 MAY 2027
S-PLEX Enhance E1	Y0550025	$\leq -10^{\circ}\text{C}$	30 NOV 2027
S-PLEX Enhance E2	Y0560024	$\leq -10^{\circ}\text{C}$	31 DEC 2027
S-PLEX Coating Reagent C1	521330	$\leq -70^{\circ}\text{C}$	30 JUN 2028
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580034	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Detect D2	Y0590026	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Enhance E3	Y0570027	$\leq -70^{\circ}\text{C}$	31 DEC 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	15.9 - 255	128	fg/mL	80% - 120%	< 20%	19.6	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		24 SEP 2025