



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IFN- β Kit

Catalog Number: K151ADRS-Series

Lot Number: K00S0104

Expiration Date: 30 APR 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IFN- β Antibody	D00T0109	2-8°C	30 APR 2026
Biotin Human IFN- β Antibody	B00T0109	2-8°C	31 MAR 2026
Human IFN- β Calibrator	A00S0068	2-8°C	31 JUL 2027
Diluent 58	M0580007	$\leq -10^{\circ}\text{C}$	30 APR 2025
S-PLEX Coating Reagent C1	351987	$\leq -70^{\circ}\text{C}$	30 APR 2024
Blocker S1	R0010811	$\leq -10^{\circ}\text{C}$	28 FEB 2025
S-PLEX Enhance E1	Y0550009	$\leq -10^{\circ}\text{C}$	28 FEB 2025
S-PLEX Enhance E2	Y0560011	$\leq -10^{\circ}\text{C}$	31 MAR 2025
S-PLEX Enhance E3	Y0570011	$\leq -70^{\circ}\text{C}$	31 MAR 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580015	$\leq -70^{\circ}\text{C}$	30 APR 2025
S-PLEX Detect D2	Y0590011	$\leq -70^{\circ}\text{C}$	31 MAR 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator	
	Recovery	CV
IFN- β	80% - 120%	< 20%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- β	16.0 - 255	255	fg/mL	80% - 120%	< 20%	32.2	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- β	1X Calibrator	196000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		12 DEC 2022