



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151ADRS-Series

Lot Number: K00S0081

Expiration Date: 28 FEB 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IFN- β Antibody	D00T0084	2-8°C	31 DEC 2025
Biotin Human IFN- β Antibody	B00T0084	2-8°C	30 SEP 2025
Human IFN- β Calibrator	A00S0031	2-8°C	31 AUG 2025
Diluent 58	M0580001	$\leq -10^{\circ}\text{C}$	28 FEB 2023
S-PLEX Coating Reagent C1	351987	$\leq -70^{\circ}\text{C}$	30 APR 2024
Blocker S1	R0010757	$\leq -10^{\circ}\text{C}$	31 JAN 2024
S-PLEX Enhance E1	Y0550008	$\leq -10^{\circ}\text{C}$	30 NOV 2024
S-PLEX Enhance E2	Y0560010	$\leq -10^{\circ}\text{C}$	31 JAN 2025
S-PLEX Enhance E3	Y0570009	$\leq -70^{\circ}\text{C}$	30 SEP 2024
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580013	$\leq -70^{\circ}\text{C}$	31 DEC 2024
S-PLEX Detect D2	Y0590009	$\leq -70^{\circ}\text{C}$	30 NOV 2024

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	63.8	fg/mL	80% - 120%	< 20%	11.7	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- β	1X Calibrator	260000	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		09 SEP 2022