



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IFN-β Kit
Catalog Number: K151ADRS-Series; K156ADSS-Series
Lot Number: K00S0131
Expiration Date: 30 Apr 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN-β Antibody	B00T0137	2-8°C	31 Aug 2026
TURBO-BOOST [™] Human IFN-β Antibody	D00T0137	2-8°C	31 Aug 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 Aug 2025
Blocker S1	R0010831	≤ -10°C	31 Jul 2025
S-PLEX Enhance E1	Y0550011	≤ -10°C	31 Jul 2025
S-PLEX Enhance E2	Y0560013	≤ -10°C	31 Aug 2025
S-PLEX Enhance E3	Y0570013	≤ -70°C	31 Jul 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580017	≤ -70°C	31 Aug 2025
S-PLEX Detect D2	Y0590013	≤ -70°C	31 Jul 2025
Human IFN-β Calibrator	A00S0081	2-8°C	29 Feb 2028
Diluent 58	M0580007	≤ -10°C	30 Apr 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	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IFN-β	7.97 - 255	15.9	fg/mL	80% - 120%	< 20%	4.77	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN-β	1X Calibrator	157000	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		27 JUL 2023