



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX® Human IFN-β Kit

**Catalog Number:** K151ADRS-Series

**Lot Number:** K00S0035

**Expiration Date:** 28 Feb 2023

#### Precision, Accuracy, and Sensitivity Specifications:

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

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IFN-β	15.9 – 255	63.8	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IFN-β	16.9	fg/mL

#### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN-β	1X Calibrator	260000	fg/mL

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IFN-β Antibody	D00T0040A	2-8°C	31 Sep 2024
Biotin Human IFN-β Antibody	B00T0040A	2-8°C	30 Apr 2024
Human IFN-β Calibrator	A00S0031	2-8°C	31 Aug 2025
Diluent 58	M0580001	≤ -10°C	28 Feb 2023
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 Aug 2023
Blocker S1	R0140004	≤ -10°C	30 Apr 2023
S-PLEX Enhance E1	Y0550005	≤ -10°C	31 May 2023
S-PLEX Enhance E2	Y0560006	≤ -10°C	31 May 2023
S-PLEX Enhance E3	Y0570005	≤ -70°C	28 Feb 2023
S-PLEX Detect D1	Y0580007	≤ -70°C	31 May 2023
S-PLEX Detect D2	Y0590005	≤ -70°C	28 Feb 2023

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		29 APR 2021