



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®]Human IFN- α 2a Kit

Catalog Number: K151P3S - Series

Lot Number: K00S0023

Expiration Date: 31 JAN 2022

Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IFN- α 2a	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IFN- α 2a	3.50 - 56.0	14.0	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IFN- α 2a	6.19	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- α 2a	750X Calibrator	44750000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IFN- α 2a Detection Antibody	D00T0027A	2-8°C	30 Nov 2023
Biotin Human IFN- α 2a Antibody	B00T0027A	2-8°C	31 Jan 2024
Human IFN- α 2a Calibrator	A00S0022	2-8°C	31 Jul 2025
Diluent 43	M0430033	\leq -10°C	31 Jan 2022
Diluent 3	M0030101	\leq -10°C	31 Dec 2022
S-PLEX Coating Reagent C1	314374	\leq -70°C	28 Feb 2023
Blocker S1	R0140003	\leq -10°C	31 Aug 2022
S-PLEX Enhance E1	Y0550003	\leq -10°C	30 Sep 2022
S-PLEX Enhance E2	Y0560005	\leq -10°C	30 Sep 2022
S-PLEX Enhance E3	Y0570003	\leq -70°C	31 Dec 2022
S-PLEX Detect D1	Y0580003	\leq -70°C	30 Nov 2022
S-PLEX Detect D2	Y0590003	\leq -70°C	31 Aug 2022

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		17 SEP 2020