



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup>Human IFN- $\alpha$ 2a Kit

**Catalog Number:** K151P3S-Series

**Lot Number:** K00S0050

**Expiration Date:** 31 MAR 2022

### Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IFN- $\alpha$ 2a	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IFN- $\alpha$ 2a	3.50 - 224	14.0	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IFN- $\alpha$ 2a	5.65	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- $\alpha$ 2a	750X Calibrator	47625000	fg/mL

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human IFN- $\alpha$ 2a Antibody	D00T0059A	2-8°C	31 OCT 2024
Biotin Human IFN- $\alpha$ 2a Antibody	B00T0059	2-8°C	30 SEP 2024
Human IFN- $\alpha$ 2a Calibrator	A00S0048	2-8°C	30 APR 2026
Diluent 43	M0430035	$\leq$ -10°C	31 MAR 2022
Diluent 3	M0030106	$\leq$ -10°C	31 AUG 2023
S-PLEX Coating Reagent C1	314374	$\leq$ -70°C	28 FEB 2023
Blocker S1	R0140004	$\leq$ -10°C	30 APR 2023
S-PLEX Enhance E1	Y0550005	$\leq$ -10°C	31 MAY 2023
S-PLEX Enhance E2	Y0560005	$\leq$ -10°C	30 SEP 2022
S-PLEX Enhance E3	Y0570004	$\leq$ -70°C	31 JAN 2023
S-PLEX Detect D1	Y0580006	$\leq$ -70°C	31 MAR 2023
S-PLEX Detect D2	Y0590005	$\leq$ -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		06 AUG 2021