



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151P3S-Series

Lot Number: K00S0070

Expiration Date: 31 AUG 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IFN- α 2a Antibody	D00T0078A	2-8°C	31 MAY 2025
Biotin Human IFN- α 2a Antibody	B00T0078A	2-8°C	31 JUL 2025
Human IFN- α 2a Calibrator	A00S0058	2-8°C	30 NOV 2026
Diluent 43	M0430047	≤ -10°C	31 MAY 2024
Diluent 3	M0030113	≤ -10°C	31 JUL 2024
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0140005	≤ -10°C	31 AUG 2023
S-PLEX Enhance E1	Y0550007	≤ -10°C	31 AUG 2024
S-PLEX Enhance E2	Y0560008	≤ -10°C	30 APR 2024
S-PLEX Enhance E3	Y0570008	≤ -70°C	30 JUN 2024
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580012	≤ -70°C	31 OCT 2024
S-PLEX Detect D2	Y0590008	≤ -70°C	31 JUL 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- α 2a	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- α 2a	3.50 - 56.0	14.0	fg/mL	80% - 120%	< 20%	5.11	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- α 2a	750X Calibrator	51000000	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		21 JUN 2022