



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151P3S-Series

Lot Number: K00S0014

Expiration Date: 31 OCT 2020

Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ				
	Range Tested	Observed	Units	Recovery	CV
IFN- α 2a	3.50 - 56.0	28.0	fg/mL	80% - 120%	< 20%

Name	LLOD	
	Observed	Units
IFN- α 2a	4.92	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- α 2a	750X Calibrator	50000000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
S-PLEX [®] 96-Well SECTOR Plate	Z00S0003	2-8°C	31 DEC 2021
TURBO-BOOST [™] Human IFN- α 2a Antibody	D00T0016A	2-8°C	31 OCT 2022
Biotin Human IFN- α 2a Antibody	B00T0016A	2-8°C	31 OCT 2022
Human IFN- α 2a Calibrator	A00S0016	2-8°C	31 JUL 2024
Diluent 43	M0430033	\leq -10°C	31 JAN 2022
Diluent 3	M0030091	\leq -10°C	31 JAN 2022
S-PLEX Coating Reagent C1	288327	\leq -70°C	31 MAY 2022
Blocker S1	R0140002	\leq -10°C	31 DEC 2021
S-PLEX Enhance E1	Y0550002	\leq -10°C	31 DEC 2021
S-PLEX Enhance E2	Y0560002	\leq -10°C	31 DEC 2021
S-PLEX Enhance E3	Y0570002	\leq -70°C	31 DEC 2021
S-PLEX Detect D1	Y0580002	\leq -70°C	31 JAN 2022
S-PLEX Detect D2	Y0590002	\leq -70°C	31 OCT 2020

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		28 JUL 2020