



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IFN- α 2a Kit
Catalog Number: K151P3S-Series
Lot Number: K00S0170
Expiration Date: 31 MAR 2026

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- α 2a Antibody	B00T0177	2-8°C	30 JUN 2027
Human IFN- α 2a Calibrator	A00S0082	2-8°C	31 MAR 2028
TURBO-BOOST™ Human IFN- α 2a Antibody	D00T0177	2-8°C	30 APR 2027
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 3	M0030126	≤ -10°C	31 JUL 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560017	≤ -10°C	31 JUL 2026
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580023	≤ -70°C	31 JUL 2026
S-PLEX Detect D2	Y0590018	≤ -70°C	31 JUL 2026
S-PLEX Enhance E3	Y0570018	≤ -70°C	30 JUN 2026

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

Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- α 2a	3.50 - 56.0	28.0	fg/mL	80% - 120%	< 20%	6.10	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- α 2a	750X Calibrator	49375000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		25 APR 2024