



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IFN- α 2a Kit
Lot Number: K00S0252
Expiration Date: 31 MAR 2026
Catalog Number: K151P3S-Series

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	A00S0048	2-8°C	30 APR 2026
TURBO-BOOST Human IFN- α 2a Antibody	D00T0177	2-8°C	30 APR 2027
S-PLEX Coating Reagent C1	446866	$\leq -70^{\circ}\text{C}$	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580029	$\leq -70^{\circ}\text{C}$	30 APR 2027
S-PLEX Detect D2	Y0590022	$\leq -70^{\circ}\text{C}$	30 APR 2027
S-PLEX Enhance E3	Y0570022	$\leq -70^{\circ}\text{C}$	30 APR 2027
Blocker S1	R0140007	$\leq -10^{\circ}\text{C}$	30 JUN 2026
Diluent 3	M0030126	$\leq -10^{\circ}\text{C}$	31 JUL 2026
Diluent 43	M0430055	$\leq -10^{\circ}\text{C}$	30 APR 2026
S-PLEX Enhance E1	Y0550020	$\leq -10^{\circ}\text{C}$	30 APR 2027
S-PLEX Enhance E2	Y0560019	$\leq -10^{\circ}\text{C}$	30 NOV 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	14.0	fg/mL	80% - 120%	< 20%	5.48	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- α 2a	750X Calibrator	47625000	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		10 FEB 2025