



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IFN- α 2a Kit
Lot Number: K00S0259
Expiration Date: 28 FEB 2027
Catalog Number: K151P3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- α 2a Antibody	B00T0244	2-8°C	31 OCT 2028
Human IFN- α 2a Calibrator	A00S0109	2-8°C	31 AUG 2029
TURBO-BOOST Human IFN- α 2a Antibody	D00T0244	2-8°C	30 SEP 2028
Blocker S1	R0140013	$\leq -10^{\circ}\text{C}$	31 JUL 2027
Diluent 3	M0030138	$\leq -10^{\circ}\text{C}$	31 AUG 2027
Diluent 43	M0430058	$\leq -10^{\circ}\text{C}$	28 FEB 2027
S-PLEX Enhance E1	Y0550023	$\leq -10^{\circ}\text{C}$	31 AUG 2027
S-PLEX Enhance E2	Y0560022	$\leq -10^{\circ}\text{C}$	31 JUL 2027
S-PLEX Coating Reagent C1	476750	$\leq -70^{\circ}\text{C}$	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580031	$\leq -70^{\circ}\text{C}$	31 JUL 2027
S-PLEX Detect D2	Y0590025	$\leq -70^{\circ}\text{C}$	30 NOV 2027
S-PLEX Enhance E3	Y0570025	$\leq -70^{\circ}\text{C}$	31 AUG 2027

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

Precision, Accuracy, and Sensitivity Specifications:

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

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- α 2a	3.50 - 56.0	14.0	fg/mL	80% - 120%	< 20%	2.92	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- α 2a	750X Calibrator	50500000	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		03 NOV 2025