



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151P3S - Series

Lot Number: K00S0136

Expiration Date: 31 Mar 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- α 2a Antibody	B00T0126	2-8°C	31 Jul 2026
TURBO-BOOST [™] Human IFN- α 2a Detection Antibody	D00T0126	2-8°C	31 Aug 2026
S-PLEX Coating Reagent C1	417723	$\leq -70^{\circ}\text{C}$	31 Aug 2025
Human IFN- α 2a Calibrator	A00S0082	2-8°C	31 Mar 2028
Diluent 43	M0430050	$\leq -10^{\circ}\text{C}$	31 Mar 2025
Diluent 3	M0030122	$\leq -10^{\circ}\text{C}$	30 Nov 2025
Blocker S1	R0010831	$\leq -10^{\circ}\text{C}$	31 Jul 2025
S-PLEX Enhance E1	Y0550011	$\leq -10^{\circ}\text{C}$	31 Jul 2025
S-PLEX Enhance E2	Y0560013	$\leq -10^{\circ}\text{C}$	31 Aug 2025
S-PLEX Enhance E3	Y0570013	$\leq -70^{\circ}\text{C}$	31 Jul 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580017	$\leq -70^{\circ}\text{C}$	31 Aug 2025
S-PLEX Detect D2	Y0590013	$\leq -70^{\circ}\text{C}$	31 Jul 2025

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	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IFN- α 2a	3.50 - 56.0	56.0	fg/mL	80% - 120%	< 20%	4.04	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- α 2a	750X Calibrator	49375000	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		02 FEB 2024