

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human IFN- α 2a Kit
Lot Number: K00S0323
Expiration Date: 28 FEB 2027
Catalog Number: K151P3S-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

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	A00S0146	2-8°C	31 JAN 2031
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™ 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:

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

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IFN- α 2a	3.50 - 56.0	7.24	fg/mL	80% - 120%	< 20%	2.84	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- α 2a	750X Calibrator	52000000	fg/mL

This lot meets MSD's non-specific binding specifications.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 APR 2026