

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human IFN- α 2a Kit
Lot Number: K00S0308
Expiration Date: 30 SEP 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	B00T0279	2-8°C	31 AUG 2029
Human IFN- α 2a Calibrator	A00S0146	2-8°C	31 JAN 2031
TURBO-BOOST Human IFN- α 2a Antibody	D00T0279	2-8°C	31 OCT 2029
Blocker S1	R0140022	$\leq -10^{\circ}\text{C}$	31 OCT 2028
Diluent 3	M0030149	$\leq -10^{\circ}\text{C}$	31 AUG 2028
Diluent 43	M0430059	$\leq -10^{\circ}\text{C}$	30 SEP 2027
S-PLEX Enhance E1	Y0550028	$\leq -10^{\circ}\text{C}$	31 MAR 2028
S-PLEX Enhance E2	Y0560027	$\leq -10^{\circ}\text{C}$	31 MAR 2028
S-PLEX Coating Reagent C1	521331	$\leq -70^{\circ}\text{C}$	31 OCT 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580037	$\leq -70^{\circ}\text{C}$	30 APR 2028
S-PLEX Detect D2	Y0590029	$\leq -70^{\circ}\text{C}$	31 MAY 2028
S-PLEX Enhance E3	Y0570030	$\leq -70^{\circ}\text{C}$	31 MAR 2028

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	19.3	fg/mL	80% - 120%	< 20%	5.67	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		22 JUN 2026