



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX[®] Human IFN- ω Kit
Lot Number: K00S0281
Expiration Date: 31 AUG 2027
Catalog Number: K151AXDS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IFN- ω Antibody	B00T0261	2-8°C	30 NOV 2028
Human IFN- ω Antibody	D00T0261	2-8°C	31 DEC 2028
Blocker S1	R0140015	$\leq -10^{\circ}\text{C}$	31 OCT 2027
S-PLEX Enhance E1	Y0550024	$\leq -10^{\circ}\text{C}$	31 OCT 2027
S-PLEX Enhance E2	Y0560023	$\leq -10^{\circ}\text{C}$	31 OCT 2027
Human IFN- ω Calibrator	A00S0136	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Coating Reagent C1	521329	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG [™] Reagent	Y0580032	$\leq -70^{\circ}\text{C}$	31 AUG 2027
S-PLEX Detect D2	Y0590026	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Enhance E3	Y0570027	$\leq -70^{\circ}\text{C}$	31 DEC 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- ω	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- ω	6.88 - 220	27.5	fg/mL	80% - 120%	< 20%	1.83	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- ω	20X Calibrator	1236000	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		20 AUG 2025