



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-12p70 Kit
Lot Number: K00S0262
Expiration Date: 31 DEC 2026
Catalog Number: K151G3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-12p70 Antibody	B00T0247	2-8°C	30 SEP 2028
Human IL-12p70 Calibrator	A00S0070	2-8°C	31 AUG 2027
TURBO-BOOST Human IL-12p70 Antibody	D00T0247	2-8°C	31 AUG 2028
Blocker S1	R0140014	≤ -10°C	30 SEP 2027
Diluent 2	M0020131	≤ -10°C	31 DEC 2026
Diluent 3	M0030138	≤ -10°C	31 AUG 2027
S-PLEX Enhance E1	Y0550022	≤ -10°C	31 JUL 2027
S-PLEX Enhance E2	Y0560022	≤ -10°C	31 JUL 2027
S-PLEX Coating Reagent C1	521329	≤ -70°C	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580031	≤ -70°C	31 JUL 2027
S-PLEX Detect D2	Y0590023	≤ -70°C	31 JUL 2027
S-PLEX Enhance E3	Y0570024	≤ -70°C	30 JUN 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	IL-12p70	80% - 120%	< 20%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-12p70	0.850 - 13.6	3.40	fg/mL	80% - 120%	< 20%	1.39	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-12p70	15X Calibrator	495000	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 JUN 2025