

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX Human IL-12p70 Kit  
**Lot Number:** K00S0206  
**Expiration Date:** 30 NOV 2025  
**Catalog Number:** K151G3S-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-12p70 Antibody	B00T0219	2-8°C	31 OCT 2027
Human IL-12p70 Calibrator	A00S0070	2-8°C	31 AUG 2027
TURBO-BOOST Human IL-12p70 Antibody	D00T0219	2-8°C	31 OCT 2027
Blocker S1	R0140009	≤ -10°C	31 OCT 2026
Diluent 2	M0020127	≤ -10°C	30 NOV 2025
Diluent 3	M0030123	≤ -10°C	31 JAN 2026
S-PLEX Enhance E1	Y0550017	≤ -10°C	31 AUG 2026
S-PLEX Enhance E2	Y0560018	≤ -10°C	31 AUG 2026
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580025	≤ -70°C	30 SEP 2026
S-PLEX Detect D2	Y0590019	≤ -70°C	30 SEP 2026
S-PLEX Enhance E3	Y0570020	≤ -70°C	31 OCT 2026

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	IL-12p70	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IL-12p70	0.850 - 13.6	13.6	fg/mL	80% - 120%	< 20%	1.99	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
<b>Review/Approval</b>	Karen Hamilla	Quality		15 JUL 2024