

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human IL-12p70 Kit
Catalog Number: K151G3S-Series
Lot Number: K00S0203
Expiration Date: 31 JAN 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-12p70 Antibody	B00T0087	2-8°C	30 NOV 2025
Human IL-12p70 Calibrator	A00S0070	2-8°C	31 AUG 2027
TURBO-BOOST™ Human IL-12p70 Antibody	D00T0087	2-8°C	31 DEC 2025
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 2	M0020123	≤ -10°C	28 FEB 2025
Diluent 3	M0030117	≤ -10°C	31 JAN 2025
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560018	≤ -10°C	31 AUG 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580021	≤ -70°C	30 APR 2026
S-PLEX Detect D2	Y0590017	≤ -70°C	30 JUN 2026
S-PLEX Enhance E3	Y0570015	≤ -70°C	30 NOV 2025

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	6.80	fg/mL	80% - 120%	< 20%	2.13	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		14 MAY 2024