



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX[®] Human IL-12p70 Kit
Lot Number: K00S0250
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	Y0550020	≤ -10°C	30 APR 2027
S-PLEX Enhance E2	Y0560020	≤ -10°C	31 MAR 2027
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG [™] Reagent	Y0580029	≤ -70°C	30 APR 2027
S-PLEX Detect D2	Y0590022	≤ -70°C	30 APR 2027
S-PLEX Enhance E3	Y0570022	≤ -70°C	30 APR 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	IL-12p70	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-12p70	0.850 - 13.6	3.40	fg/mL	80% - 120%	< 20%	1.60	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.

Review/Approval	Karen Hamilla	Quality		31 DEC 2024
------------------------	---------------	---------	---	-------------