



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL12p70 Kit

Catalog Number: K151G3S-Series

Lot Number: K00S0117

Expiration Date: 30 APR 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IL12p70 Antibody	D00T0087	2-8°C	31 DEC 2025
Biotin Human IL12p70 Antibody	B00T0087	2-8°C	30 NOV 2025
Human IL-12p70 Calibrator	A00S0070	2-8°C	31 AUG 2027
Diluent 2	M0020123	≤ -10°C	28 FEB 2025
Diluent 3	M0030117	≤ -10°C	31 JAN 2025
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0010768	≤ -10°C	31 MAY 2024
S-PLEX Enhance E1	Y0550009	≤ -10°C	28 FEB 2025
S-PLEX Enhance E2	Y0560010	≤ -10°C	31 JAN 2025
S-PLEX Enhance E3	Y0570011	≤ -70°C	31 MAR 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580014	≤ -70°C	28 FEB 2025
S-PLEX Detect D2	Y0590009	≤ -70°C	30 NOV 2024

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator	
	Recovery	CV
IL-12p70	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-12p70	0.850 - 13.6	13.6	fg/mL	80% - 120%	< 20%	1.61	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-12p70	15X Calibrator	495000	fg/mL

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 DEC 2022