



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-12p70 Kit

**Catalog Number:** K151G3S-Series

**Lot Number:** K00S0060

**Expiration Date:** 31 AUG 2023

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL12p70 Antibody	D00T0071	2-8°C	31 DEC 2024
Biotin Human IL12p70 Antibody	B00T0071	2-8°C	30 NOV 2024
Human IL-12p70 Calibrator	A00S0044	2-8°C	31 DEC 2025
Diluent 2	M0020117	≤ -10°C	31 OCT 2023
Diluent 3	M0030109	≤ -10°C	31 JAN 2024
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 Aug 2023
Blocker S1	R0140005	≤ -10°C	31 Aug 2023
S-PLEX Enhance E1	R0140006	≤ -10°C	31 Jan 2024
S-PLEX Enhance E2	Y0560007	≤ -10°C	31 Oct 2023
S-PLEX Enhance E3	Y0570007	≤ -70°C	29 Feb 2024
S-PLEX Detect D1	Y0580009	≤ -70°C	29 Feb 2024
S-PLEX Detect D2	Y0590007	≤ -70°C	29 Feb 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	6.80	fg/mL	80% - 120%	< 20%	1.47	fg/mL

#### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-12p70	15X Calibrator	471000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		27 JAN 2022