



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX® Human IL-12p70 Kit

**Catalog Number:** K151G3S-Series

**Lot Number:** K00S0043

**Expiration Date:** 28 FEB 2022

#### Precision, Accuracy, and Sensitivity Specifications:

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

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

LLOD		
Name	Observed	Units
IL-12p70	1.85	fg/mL

#### Stock Concentration of Calibrator:

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

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-12p70 Antibody	D00T0047A	2-8°C	31 Jul 2024
Biotin Human IL-12p70 Antibody	B00T0047A	2-8°C	31 May 2024
Human IL-12p70 Calibrator	A00S0044	2-8°C	31 Dec 2025
Diluent 2	M0020108	≤ -10°C	28 Feb 2022
Diluent 3	M0030104	≤ -10°C	30 Jun 2023
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 Aug 2023
Blocker S1	R0140004	≤ -10°C	30 Apr 2023
S-PLEX Enhance E1	Y0550004	≤ -10°C	28 Feb 2023
S-PLEX Enhance E2	Y0560006	≤ -10°C	31 May 2023
S-PLEX Enhance E3	Y0570004	≤ -70°C	31 Jan 2023
S-PLEX Detect D1	Y0580006	≤ -70°C	31 Mar 2023
S-PLEX Detect D2	Y0590004	≤ -70°C	31 Jan 2023

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 MAR 2021