



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-5 Kit  
**Catalog Number:** K151J3S-Series; K156J3S-Series  
**Lot Number:** K00S0026  
**Expiration Date:** 31 Jan 2022

### Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IL-5	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-5	3.00 – 48.0	48.0	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IL-5	1.91	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-5	30X Calibrator	962000	fg/mL

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-5 Antibody	D00T0030A	2-8°C	30 Nov 2023
Biotin Human IL-5 Antibody	B00T0030A	2-8°C	31 Mar 2024
Human IL-5 Calibrator	A00S0008	2-8°C	31 Jul 2024
Diluent 43	M0430033	≤ -10°C	31 Jan 2022
Diluent 3	M0030104	≤ -10°C	30 Jun 2023
S-PLEX Coating Reagent C1	314374	≤ -70°C	28 Feb 2023
Blocker S1	R0140004	≤ -10°C	30 Apr 2023
S-PLEX Enhance E1	Y0550004	≤ -10°C	28 Feb 2023
S-PLEX Enhance E2	Y0560004	≤ -10°C	31 Jan 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		04 FEB 2021