



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-4 Kit

**Catalog Number:** K151A3S-Series

**Lot Number:** K00S0025

**Expiration Date:** 31 JAN 2023

### Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IL-4	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-4	1.23 - 19.6	9.80	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IL-4	0.552	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-4	20X Calibrator	179000	fg/mL

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-4 Antibody	D00T0029A	2-8°C	30 Apr 2024
Biotin Human IL-4 Antibody	B00T0029A	2-8°C	31 Jan 2024
Human IL-4 Calibrator	A00S0020	2-8°C	30 Sep 2024
Diluent 2	M0020114	≤ -10°C	31 Mar 2023
Diluent 3	M0030103	≤ -10°C	30 Apr 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	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
<b>Review/Approval</b>	Karen Hamilla	Quality		29 JAN 2021