



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX® Human IL-17A Kit

**Catalog Number:** K151C3S-Series

**Lot Number:** K00S0044

**Expiration Date:** 28 FEB 2023

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

Calibrator		
Name	Recovery	CV
IL-17A	75% - 125%	< 25%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-17A	3.75 - 120	120	fg/mL	75% - 125%	< 25%

LLOD		
Name	Observed	Units
IL-17A	12.0	fg/mL

#### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-17A	10X Calibrator	2660000	fg/mL

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-17A Antibody	D00T0048A	2-8°C	30 SEP 2024
Biotin Human IL-17A Antibody	B00T0048A	2-8°C	30 APR 2024
Human IL-17A Calibrator	A00S0023	2-8°C	31 Jul 2025
Diluent 43	M0430044	≤ -10°C	30 Sep 2023
Diluent 3	M0030106	≤ -10°C	31 Aug 2023
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 Aug 2023
Blocker S1	R0140004	≤ -10°C	30 Apr 2023
S-PLEX Enhance E1	Y0550005	≤ -10°C	31 May 2023
S-PLEX Enhance E2	Y0560006	≤ -10°C	31 May 2023
S-PLEX Enhance E3	Y0570005	≤ -70°C	28 Feb 2023
S-PLEX Detect D1	Y0580007	≤ -70°C	31 May 2023
S-PLEX Detect D2	Y0590005	≤ -70°C	28 Feb 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		27 MAY 2021