



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-17A Kit

**Catalog Number:** K151C3S-Series

**Lot Number:** K00S0011A

**Expiration Date:** 30 APR 2021

### Precision, Accuracy, and Sensitivity Specifications:

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

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-17A	3.75-60.0	30.0	fg/mL	75% - 125%	< 25%

LLOD		
Name	Observed	Units
IL-17A	13.36	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-17A	10X Calibrator	2350000	fg/mL

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-17A Antibody	D00T0005A	2-8°C	31 OCT 2022
Biotin Human IL-17A Antibody	B00T0005A	2-8°C	31 OCT 2022
Human IL-17A Calibrator	A00S0018	2-8°C	30 SEP 2024
Diluent 43	M0430033	≤ -10°C	31 JAN 2022
Diluent 3	M0030091	≤ -10°C	31 JAN 2022
S-PLEX Coating Reagent C1	265437	≤ -70°C	30 SEP 2021
Blocker S1	R0140001	≤ -10°C	31 MAY 2021
S-PLEX Enhance E1	Y0550002	≤ -10°C	31 DEC 2021
S-PLEX Enhance E2	Y0560002	≤ -10°C	31 DEC 2021
S-PLEX Enhance E3	Y0570001	≤ -70°C	31 MAY 2021
S-PLEX Detect D1	Y0580002	≤ -70°C	31 JAN 2022
S-PLEX Detect D2	Y0590001	≤ -70°C	30 APR 2021

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		22 FEB 2021