



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-17A Kit
Catalog Number: K151C3S-Series; K156C3S-Series
Lot Number: K00S0016
Expiration Date: 31 Jan 2022

Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
Analyte	75% - 125%	< 25%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-17A	3.75 - 240	120	fg/mL	75% - 125%	< 25%

LLOD		
Name	Observed	Units
IL-17A	57.8	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	D00T0024A	2-8°C	30 Sep 2023
Biotin Human IL-17A Antibody	B00T0024A	2-8°C	31 Dec 2023
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	314374	≤ -70°C	28 Feb 2023
Blocker S1	R0140003	≤ -10°C	31 Aug 2022
S-PLEX Enhance E1	Y0550003	≤ -10°C	30 Sep 2022
S-PLEX Enhance E2	Y0560003	≤ -10°C	31 Mar 2022
S-PLEX Enhance E3	Y0570003	≤ -70°C	31 Dec 2022
S-PLEX Detect D1	Y0580003	≤ -70°C	30 Nov 2022
S-PLEX Detect D2	Y0590003	≤ -70°C	31 Aug 2022

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 OCT 2020