



EXTENSION MEMO

Page 1 of 1

Product Description(s): S-PLEX® Human IL-17A Kit


Catalog Number(s) / Size(s): K151C3S-Series; K156C3S-Series

Lot Number: K00S0071

Based on additional testing, the original date of expiration for this kit has been extended by 11 months. The expiration date of K00S0071 has been extended to 31 JUL 2024.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 43	M0430047	31 MAY 2024	31 JUL 2024
S-PLEX Coating Reagent C1	351987	30 APR 2024	31 JUL 2024
Blocker S1	R0140005	31 AUG 2023	31 JUL 2024
S-PLEX Enhance E2	Y0560008	30 APR 2024	31 JUL 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		19 JUL 2023

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX® Human IL-17A Kit
Catalog Number: K151C3S-Series; K156C3S-Series
Lot Number: K00S0071
Expiration Date: 31 AUG 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-17A Antibody	D00T0079	2-8°C	31 MAY 2025
Biotin Human IL-17A Antibody	B00T0079	2-8°C	30 SEP 2025
Human IL-17A Calibrator	A00S0045	2-8°C	31 DEC 2025
Diluent 43	M0430047	≤ -10°C	31 MAY 2024
Diluent 3	M0030113	≤ -10°C	31 JUL 2024
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0140005	≤ -10°C	31 AUG 2023
S-PLEX Enhance E1	Y0550007	≤ -10°C	31 AUG 2024
S-PLEX Enhance E2	Y0560008	≤ -10°C	30 APR 2024
S-PLEX Enhance E3	Y0570009	≤ -70°C	30 SEP 2024
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580012	≤ -70°C	31 OCT 2024
S-PLEX Detect D2	Y0590008	≤ -70°C	31 JUL 2024

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-17A	7.50 - 120	60.0	fg/mL	75% - 125%	< 25%	11.3	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-17A	10X Calibrator	2400000	fg/mL

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 JUL 2022