



## EXTENSION MEMO

Page 1 of 1

**Product Description(s):** S-PLEX® Human IL-21 Kit

**Catalog Number(s) / Size(s):** K151ADNS-Series

**Lot Number:** K00S0127

Based on additional testing, the original date of expiration for this Kit has been extended by 12 months. The expiration date of K00S0127 has been extended to 28 FEB 2025.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
TURBO-BOOST™ Human IL-21 Antibody	D00T0042A	31 MAY 2024	28 FEB 2025
Biotin Human IL-21 Antibody	B00T0042A	30 APR 2024	28 FEB 2025
S-PLEX Coating Reagent C1	351987	30 APR 2024	28 FEB 2025
Blocker S1	R0010757	29 FEB 2024	28 FEB 2025
S-PLEX Enhance E1	Y0550007	31 AUG 2024	28 FEB 2025
S-PLEX Enhance E2	Y0560008	30 APR 2024	28 FEB 2025
S-PLEX Enhance E3	Y0570009	30 SEP 2024	28 FEB 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580012	31 OCT 2024	28 FEB 2025
S-PLEX Detect D2	Y0590008	31 JUL 2024	28 FEB 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		12 JAN 2024

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



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-21 Kit

**Catalog Number:** K151ADNS-Series

**Lot Number:** K00S0127

**Expiration Date:** 29 FEB 2024

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human IL-21 Antibody	D00T0042A	2-8°C	31 MAY 2024
Biotin Human IL-21 Antibody	B00T0042A	2-8°C	30 APR 2024
Human IL-21 Calibrator	A00S0072	2-8°C	31 AUG 2027
Diluent 2	M0020124	≤ -10°C	28 FEB 2025
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0010757	≤ -10°C	29 FEB 2024
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 <sup>™</sup> 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-21	80% - 120%	< 20%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-21	18.8 - 300	18.8	fg/mL	80% - 120%	< 20%	9.91	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-21	1X Calibrator	86700	fg/mL

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		03 MAR 2023