



**RFE MEMO**

Page 1 of 1

**Product Description(s):** S-PLEX® Human IL-22 Kit

**Catalog Number(s) / Size(s):** K151H3S - Series

**Lot Number:** K00S0018

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0018 has been extended to 31 JAN 2024.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
TURBO-BOOST™ Human IL-22 Antibody	D00T0025A	30 SEP 2023	31 JAN 2024
Biotin Human IL-22 Antibody	B00T0025A	30 NOV 2023	31 JAN 2024
Diluent 43	M0430044	30 SEP 2023	31 JAN 2024
Diluent 3	M0030106	31 AUG 2023	31 JAN 2024
S-PLEX Coating Reagent C1	329998	31 AUG 2023	31 JAN 2024
Blocker S1	R0140004	30 APR 2023	31 JAN 2024
S-PLEX Enhance E1	Y0550005	31 MAY 2023	31 JAN 2024
S-PLEX Enhance E2	Y0560006	31 MAY 2023	31 JAN 2024
S-PLEX Enhance E3	Y0570004	31 JAN 2023	31 JAN 2024
S-PLEX Detect D1	Y0580006	31 MAR 2023	31 JAN 2024
S-PLEX Detect D2	Y0590004	31 JAN 2023	31 JAN 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		16 DEC 2022

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-22 Kit

**Catalog Number:** K151H3S-Series

**Lot Number:** K00S0018

**Expiration Date:** 31 JAN 2023

### Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IL-22	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-22	3.75 - 60.0	15.0	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IL-22	2.67	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-22	15X Calibrator	537000	fg/mL

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human IL-22 Antibody	D00T0025A	2-8°C	30 Sep 2023
Biotin Human IL-22 Antibody	B00T0025A	2-8°C	30 Nov 2023
Human IL-22 Calibrator	A00S0025	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	Y0570004	≤ -70°C	31 Jan 2023
S-PLEX Detect D1	Y0580006	≤ -70°C	31 Mar 2023
S-PLEX Detect D2	Y0590004	≤ -70°C	31 Jan 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		18 NOV 2021