



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** K151H3S - Series

**Lot Number:** K00S0020

**Expiration Date:** 31 MAY 2021

#### 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	7.50	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IL-22	2.11	fg/mL

#### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-22	15X Calibrator	532000	fg/mL

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-22 Antibody	D00T0010A	2-8°C	31 Oct 2022
Biotin Human IL-22 Antibody	B00T0010A	2-8°C	31 Oct 2022
Human IL-22 Calibrator	A00S0014	2-8°C	31 Jul 2024
Diluent 43	M0430033	≤ -10°C	31 Jan 2022
Diluent 3	M0030091	≤ -10°C	31 Jan 2022
S-PLEX Coating Reagent C1	288327	≤ -70°C	31 May 2022
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	Y0570002	≤ -70°C	31 Dec 2021
S-PLEX Detect D1	Y0580002	≤ -70°C	31 Jan 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		14 AUG 2020