



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®]Human IL-2 Kit
Catalog Number: K151Z2S-Series; K156Z2S-Series
Lot Number: K00S0024
Expiration Date: 31 AUG 2022

Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IL-2	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-2	2.75 - 44.0	44.0	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IL-2	6.48	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-2	30X Calibrator	1420000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-2 Antibody	D00T0028A	2-8°C	30 Nov 2023
Biotin Human IL-2 Antibody	B00T0028A	2-8°C	31 Mar 2024
Human IL-2 Calibrator	A00S0001	2-8°C	31 Jul 2024
Diluent 2	M0020113	≤ -10°C	30 Nov 2022
Diluent 3	M0030102	≤ -10°C	31 Jan 2023
S-PLEX Coating Reagent C1	314374	≤ -70°C	28 Feb 2023
Blocker S1	R0140003	≤ -10°C	31 Aug 2022
S-PLEX Enhance E1	Y0550004	≤ -10°C	28 Feb 2023
S-PLEX Enhance E2	Y0560004	≤ -10°C	31 Jan 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	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		12 NOV 2020