



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-2 Kit  
**Catalog Number:** K151Z2S-Series; K156Z2S-Series  
**Lot Number:** K00S0001A  
**Expiration Date:** 30 APR 2021

#### Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ				
	Range Tested	Observed	Units	Recovery	CV
IL-2	2.75 – 44.0	22.0	fg/mL	80% - 120%	< 20%

Name	LLOD	
	Observed	Units
IL-2	7.32	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	D00T0002A	2-8°C	31 OCT 2022
Biotin Human IL-2 Antibody	B00T0002A	2-8°C	31 OCT 2022
Human IL-2 Calibrator	A00S0001	2-8°C	31 JUL 2024
Diluent 2	M0020107	≤ -10°C	31 JAN 2022
Diluent 3	M0030091	≤ -10°C	31 JAN 2022
S-PLEX Coating Reagent C1	265437	≤ -70°C	30 SEP 2021
Blocker S1	R0140001	≤ -10°C	31 MAY 2021
S-PLEX Enhance E1	Y0550001	≤ -10°C	30 APR 2021
S-PLEX Enhance E2	Y0560001	≤ -10°C	30 APR 2021
S-PLEX Enhance E3	Y0570001	≤ -70°C	31 MAY 2021
S-PLEX Detect D1	Y0580001	≤ -70°C	30 APR 2021
S-PLEX Detect D2	Y0590001	≤ -70°C	30 APR 2021

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		28 JUL 2020