



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-10 Kit

Catalog Number: K151Y2S - Series

Lot Number: K00S0015

Expiration Date: 31 DEC 2021

Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IL-10	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-10	2.45 - 39.2	4.9	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IL-10	1.56	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-10	20X Calibrator	380000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IL-10 Antibody	D00T0022A	2-8°C	30 SEP 2023
Biotin Human IL-10 Antibody	B00T0022A	2-8°C	31 OCT 2023
Human IL-10 Calibrator	A00S0004	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	314374	≤ -70°C	28 FEB 2023
Blocker S1	R0140002	≤ -10°C	31 DEC 2021
S-PLEX Enhance E1	Y0550003	≤ -10°C	30 SEP 2022
S-PLEX Enhance E2	Y0560003	≤ -10°C	31 MAR 2022
S-PLEX Enhance E3	Y0570002	≤ -70°C	31 DEC 2021
S-PLEX Detect D1	Y0580003	≤ -70°C	30 NOV 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