



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151Y2S-Series

Lot Number: K00S0049

Expiration Date: 31 Jul 2022

Precision, Accuracy, and Sensitivity Specifications:

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

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

Name	LLOD	
	Observed	Units
IL-10	0.970	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	D00T0046A	2-8°C	31 Jul 2024
Biotin Human IL-10 Antibody	B00T0046A	2-8°C	30 Apr 2024
Human IL-10 Calibrator	A00S0004	2-8°C	31 Jul 2024
Diluent 2	M0020112	≤ -10°C	31 Jul 2022
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	Y0570005	≤ -70°C	28 Feb 2023
S-PLEX Detect D1	Y0580007	≤ -70°C	31 May 2023
S-PLEX Detect D2	Y0590005	≤ -70°C	28 Feb 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
Review/Approval	Karen Hamilla	Quality		16 NOV 2021