



RFE MEMO

Page 1 of 1

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

Catalog Number(s) / Size(s): K151Y2S-Series

Lot Number: K00S0078

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0078 has been extended to 31 JUL 2023.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 2	M0020112	31 JUL 2022	31 JUL 2023
Blocker S1	R0140004	30 APR 2023	31 JUL 2023
S-PLEX Enhance E1	Y0550005	31 May 2023	31 JUL 2023
S-PLEX Enhance E2	Y0560006	31 May 2023	31 JUL 2023
S-PLEX Enhance E3	Y0570005	28 Feb 2023	31 JUL 2023
S-PLEX Detect D2	Y0590005	28 Feb 2023	31 JUL 2023

Function	Name	Signature	Date
Quality	Karen Hamilla		13 APR 2022

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151Y2S-Series

Lot Number: K00S0078

Expiration Date: 31 JUL 2022

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	Y0580011	≤ -70°C	31 May 2024
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.

Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-10	2.45 - 39.2	4.90	fg/mL	80% - 120%	< 20%	1.20	fg/mL

Stock Concentration of Calibrator:

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

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		13 APR 2022