



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-10 Kit

**Catalog Number:** K151Y2S-Series

**Lot Number:** K00S0083

**Expiration Date:** 31 JAN 2024

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human IL-10 Antibody	D00T0086	2-8°C	30 NOV 2025
Biotin Human IL-10 Antibody	B00T0086	2-8°C	30 SEP 2025
Human IL-10 Calibrator	A00S0066	2-8°C	31 AUG 2027
Diluent 2	M0020122	≤ -10°C	31 OCT 2024
Diluent 3	M0030114	≤ -10°C	30 SEP 2024
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0010757	≤ -10°C	31 JAN 2024
S-PLEX Enhance E1	Y0550008	≤ -10°C	30 NOV 2024
S-PLEX Enhance E2	Y0560010	≤ -10°C	31 JAN 2025
S-PLEX Enhance E3	Y0570009	≤ -70°C	30 SEP 2024
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580013	≤ -70°C	31 DEC 2024
S-PLEX Detect D2	Y0590009	≤ -70°C	30 NOV 2024

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	9.80	fg/mL	80% - 120%	< 20%	2.28	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-10	20X Calibrator	398000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		10 NOV 2022