



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-6 Kit

**Catalog Number:** K151B3S-Series

**Lot Number:** K00S0086

**Expiration Date:** 30 APR 2024

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human IL-6 Antibody	D00T0089	2-8°C	30 SEP 2025
Biotin Human IL-6 Antibody	B00T0089	2-8°C	31 JAN 2026
Human IL-6 Calibrator	A00S0021	2-8°C	30 JUN 2025
Diluent 2	M0020123	≤ -10°C	28 FEB 2025
Diluent 3	M0030117	≤ -10°C	31 JAN 2025
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0010768	≤ -10°C	31 MAY 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	Y0570010	≤ -70°C	31 JAN 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580013	≤ -70°C	31 DEC 2024
S-PLEX Detect D2	Y0590010	≤ -70°C	28 FEB 2025

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-6	80% - 120%	< 20%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-6	0.325 – 5.20	5.20	fg/mL	80% - 120%	< 20%	1.25	fg/mL

#### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-6	120X Calibrator	768000	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		28 OCT 2022