



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-6 Kit
Catalog Number: K151B3S-Series; K156B3S-Series
Lot Number: K00S0065
Expiration Date: 31 AUG 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IL-6 Detection Antibody	D00T0074	2-8°C	31 JUL 2025
Biotin Human IL-6 Capture Antibody	B00T0074	2-8°C	31 JAN 2025
Human IL-6 Calibrator	A00S0021	2-8°C	30 JUN 2025
Diluent 2	M0020117	≤ -10°C	31 OCT 2023
Diluent 3	M0030111	≤ -10°C	31 MAY 2024
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 AUG 2023
Blocker S1	R0140005	≤ -10°C	31 AUG 2023
S-PLEX Enhance E1	R0140006	≤ -10°C	31 JAN 2024
S-PLEX Enhance E2	Y0560007	≤ -10°C	31 OCT 2023
S-PLEX Enhance E3	Y0570007	≤ -70°C	29 FEB 2024
S-PLEX Detect D1	Y0580009	≤ -70°C	29 FEB 2024
S-PLEX Detect D2	Y0590007	≤ -70°C	29 FEB 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-6	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IL-6	0.325 - 5.20	5.20	fg/mL	80% - 120%	< 20%	0.718	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
Review/Approval	Karen Hamilla	Quality		08 MAR 2022