



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®]Human IL-3 Kit

Catalog Number: K151ADQS - Series

Lot Number: K00S0169

Expiration Date: 30 APR 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-3 Antibody	B00T0043A	2-8°C	30 APR 2024
TURBO-BOOST [™] Human IL-3 Antibody	D00T0043A	2-8°C	31 MAY 2024
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
Blocker S1	R0010811	≤ -10°C	28 FEB 2025
S-PLEX Enhance E1	Y0550013	≤ -10°C	30 NOV 2025
S-PLEX Enhance E2	Y0560011	≤ -10°C	31 MAR 2025
S-PLEX Enhance E3	Y0570010	≤ -70°C	31 JAN 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580015	≤ -70°C	30 APR 2025
S-PLEX Detect D2	Y0590011	≤ -70°C	31 MAR 2025
Human IL-3 Calibrator	A00S0033	2-8°C	31 AUG 2025
Diluent 58	M0580009	≤ -10°C	30 NOV 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-3	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IL-3	10.6 - 170	42.5	fg/mL	80% - 120%	< 20%	4.11	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-3	1X Calibrator	38000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		21 NOV 2023