



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-5 Kit
Catalog Number: K151J3S-Series; K156J3S-Series
Lot Number: K00S0008A
Expiration Date: 30 APR 2021

Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator	
	Recovery	CV
IL-5	80% - 120%	< 20%

Name	LLOQ				
	Range Tested	Observed	Units	Recovery	CV
IL-5	3.00 - 48.0	6.00	fg/mL	80% - 120%	< 20%

Name	LLOD	
	Observed	Units
IL-5	2.02	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-5	30X Calibrator	962000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IL-5 Antibody	D00T0011A	2-8°C	31 OCT 2022
Biotin Human IL-5 Antibody	B00T0011A	2-8°C	31 OCT 2022
Human IL-5 Calibrator	A00S0008	2-8°C	31 JUL 2024
Diluent 43	M0430032	≤ -10°C	30 SEP 2021
Diluent 3	M0030091	≤ -10°C	31 JAN 2022
S-PLEX Coating Reagent C1	265437	≤ -70°C	30 SEP 2021
Blocker S1	R0140001	≤ -10°C	31 MAY 2021
S-PLEX Enhance E1	Y0550001	≤ -10°C	30 APR 2021
S-PLEX Enhance E2	Y0560002	≤ -10°C	31 DEC 2021
S-PLEX Enhance E3	Y0570001	≤ -70°C	31 MAY 2021
S-PLEX Detect D1	Y0580001	≤ -70°C	30 APR 2021
S-PLEX Detect D2	Y0590001	≤ -70°C	30 APR 2021

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 JUL 2020