



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-9 Kit

Catalog Number: K151R3S - Series

Lot Number: K00S0009A

Expiration Date: 30 APR 2021

Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IL-9	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-9	9.75 - 156	19.5	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IL-9	8.57	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-9	15X Calibrator	1600000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-9 Antibody	D00T0018A	2-8°C	31 Oct 2022
Biotin Human IL-9 Antibody	B00T0018A	2-8°C	31 Oct 2022
Human IL-9 Calibrator	A00S0010	2-8°C	31 Jul 2024
Diluent 43	M0430033	≤ -10°C	31 Jan 2022
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		19 MAY 2020