



RFE MEMO

Page 1 of 1

Product Description(s): S-PLEX® Human IL-9 Kit

Catalog Number(s) / Size(s): K151R3S-Series

Lot Number: K00S0089

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0089 has been extended to 31 OCT 2023.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
TURBO-BOOST™ Human IL-9 Antibody	D00T0018A	31 OCT 2022	31 OCT 2023
Biotin Human IL-9 Antibody	B00T0018A	31 OCT 2022	31 OCT 2023
S-PLEX Detect D1	Y0580006	31 MAR 2023	31 OCT 2023

Function	Name	Signature	Date
Quality	Karen Hamilla		21 JUN 2022

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151R3S-Series

Lot Number: K00S0089

Expiration Date: 31 OCT 2022

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	M0430047	≤ -10°C	31 MAY 2024
Diluent 3	M0030113	≤ -10°C	31 JUL 2024
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R010757	≤ -10°C	31 JAN 2024
S-PLEX Enhance E1	Y0550007	≤ -10°C	31 AUG 2024
S-PLEX Enhance E2	Y0560007	≤ -10°C	31 OCT 2023
S-PLEX Enhance E3	Y0570008	≤ -70°C	30 JUN 2024
S-PLEX Detect D1	Y0580006	≤ -70°C	31 MAR 2023
S-PLEX Detect D2	Y0590006	≤ -70°C	31 OCT 2023

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-9	75% - 125%	< 25%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-9	9.75 - 156	19.5	fg/mL	75% - 125%	< 25%	8.03	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-9	15X Calibrator	1600000	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		21 JUN 2022