



## EXTENSION MEMO

Page 1 of 1

**Product Description(s):** S-PLEX® SARS-CoV-2 N Kit


**Catalog Number(s) / Size(s):** K150ADHS-Series

**Lot Number:** K00S0188

Based on additional testing, the original date of expiration for this Kit has been extended by 24 months. The expiration date of K00S0188 has been extended to 31 MAR 2025.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Biotin SARS-CoV-2 N Antibody	B00T0045A	31 MAR 2024	31 MAR 2025
TURBO-BOOST™ SARS-CoV-2 N Antibody	D00T0045A	31 MAR 2024	31 MAR 2025
SARS-CoV-2 N Calibrator	A00S0034	31 MAR 2023	31 MAR 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		13 MAR 2024

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

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> SARS-CoV-2 N Kit  
**Catalog Number:** K150ADHS-Series  
**Lot Number:** K00S0188  
**Expiration Date:** 31 MAR 2023

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin SARS-CoV-2 N Antibody	B00T0045A	2-8°C	31 MAR 2024
TURBO-BOOST <sup>™</sup> SARS-CoV-2 N Antibody	D00T0045A	2-8°C	31 MAR 2024
Blocker D-R	R0060030	≤ -10°C	31 MAR 2026
Blocker S1	R0010831	≤ -10°C	31 JUL 2025
Diluent 6l	M0610004	≤ -10°C	31 AUG 2025
S-PLEX Enhance E1	Y0550012	≤ -10°C	31 AUG 2025
S-PLEX Enhance E2	Y0560013	≤ -10°C	31 AUG 2025
SARS-CoV-2 N Calibrator	A00S0034	≤ -70°C	31 MAR 2023
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580017	≤ -70°C	31 AUG 2025
S-PLEX Detect D2	Y0590013	≤ -70°C	31 JUL 2025
S-PLEX Enhance E3	Y0570013	≤ -70°C	31 JUL 2025

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

**Precision, Accuracy, and Sensitivity Specifications:**

		Calibrator	
Spot Number	Name	Recovery	CV
1	SARS-CoV-2 N	75% - 125%	< 25%

		LLOQ				LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
SARS-CoV-2 N	86.3 - 1380	173	fg/mL	75% - 125%	< 25%	57.8	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
SARS-CoV-2 N	20X Calibrator	21800000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		13 MAR 2024