



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX[®] SARS-CoV-2 N Kit
Lot Number: K00S0189
Expiration Date: 31 AUG 2025
Catalog Number: K150ADHS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin SARS-CoV-2 N Antibody	B00T0198	2-8°C	31 DEC 2027
TURBO-BOOST SARS-CoV-2 N Antibody	D00T0225	2-8°C	31 JAN 2028
Blocker S1	R0140010	≤ -10°C	31 DEC 2026
Diluent 61	M0610004	≤ -10°C	31 AUG 2025
S-PLEX Enhance E1	Y0550018	≤ -10°C	30 NOV 2026
S-PLEX Enhance E2	Y0560019	≤ -10°C	30 NOV 2026
SARS-CoV-2 N Calibrator	A00S0104	≤ -70°C	30 SEP 2026
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG [™] Reagent	Y0580027	≤ -70°C	31 OCT 2026
S-PLEX Detect D2	Y0590020	≤ -70°C	30 NOV 2026
S-PLEX Enhance E3	Y0570020	≤ -70°C	31 OCT 2026

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

Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
SARS CoV-2 N	86.3 - 1380	690	fg/mL	75% - 125%	< 25%	92.4	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
SARS CoV-2 N	20X Calibrator	23200000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 NOV 2024