



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX[®] SARS-CoV-2 N Kit
Lot Number: K00S0240
Expiration Date: 30 SEP 2026
Catalog Number: K150ADHS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin SARS-CoV-2 N Antibody	B00T0240	2-8°C	30 APR 2028
TURBO-BOOST SARS-CoV-2 N Antibody	D00T0240	2-8°C	31 AUG 2028
Blocker S1	R0140013	≤ -10°C	31 JUL 2027
Diluent 61	M0610005	≤ -10°C	30 NOV 2027
S-PLEX Enhance E1	Y0550022	≤ -10°C	31 JUL 2027
S-PLEX Enhance E2	Y0560021	≤ -10°C	30 APR 2027
SARS-CoV-2 N Calibrator	A00S0104	≤ -70°C	30 SEP 2026
S-PLEX Coating Reagent C1	521329	≤ -70°C	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG [™] Reagent	Y0580030	≤ -70°C	30 APR 2027
S-PLEX Detect D2	Y0590023	≤ -70°C	31 JUL 2027
S-PLEX Enhance E3	Y0570024	≤ -70°C	30 JUN 2027

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	345	fg/mL	75% - 125%	< 25%	88.6	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		18 JUL 2025