



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX® SARS-CoV-2 N Kit

Catalog Number: K150ADHS-Series

Lot Number: K00S0031

Expiration Date: 31 AUG 2022

Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ				
	Range Tested	Observed	Units	Recovery	CV
SARS-CoV-2 N	86.3-1380	172.5	fg/mL	75% - 125%	< 25%

Name	LLOD	
	Observed	Units
SARS-CoV-2 N	101	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
SARS-CoV-2 N	20X Calibrator	19220000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ SARS-CoV-2 N Antibody	D00T0035A	2-8°C	31 JAN 2024
Biotin SARS-CoV-2 N Antibody	B00T0035A	2-8°C	31 JAN 2024
SARS-CoV-2 N Calibrator	A00V0002	≤ -70°C	31 JAN 2023
S-PLEX Coating Reagent C1	314374	≤ -70°C	28 FEB 2023
Blocker S1	R0140003	≤ -10°C	31 AUG 2022
Blocker D-R	R0060024	≤ -10°C	31 DEC 2022
S-PLEX Enhance E1	Y0550003	≤ -10°C	30 SEP 2022
S-PLEX Enhance E2	Y0560005	≤ -10°C	30 SEP 2022
S-PLEX Enhance E3	Y0570003	≤ -70°C	31 DEC 2022
S-PLEX Detect D1	Y0580003	≤ -70°C	30 NOV 2022
S-PLEX Detect D2	Y0590003	≤ -70°C	31 AUG 2022

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		17 SEP 2020