



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K150ADHS-Series

Lot Number: K00S0038

Expiration Date: 31 DEC 2021

Precision, Accuracy, and Sensitivity Specifications:

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

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

LLOD		
Name	Observed	Units
SARS-CoV-2 N	52.5	fg/mL

Stock Concentration of Calibrator:

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

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] SARS-CoV-2 N Antibody	D00T0045A	2-8°C	31 Mar 2024
Biotin SARS-CoV-2 N Antibody	B00T0045A	2-8°C	31 Mar 2024
SARS-CoV-2 N Calibrator	A00S0034	≤ -70°C	31 Mar 2023
Diluent 61	M0610001	≤ -10°C	31 Mar 2023
S-PLEX Coating Reagent C1	288327	≤ -70°C	31 May 2022
Blocker S1	R0140002	≤ -10°C	31 Dec 2021
S-PLEX Enhance E1	Y0550004	≤ -10°C	28 Feb 2023
S-PLEX Enhance E2	Y0560004	≤ -10°C	31 Jan 2023
S-PLEX Enhance E3	Y0570003	≤ -70°C	31 Dec 2022
S-PLEX Detect D1	Y0580005	≤ -70°C	31 Mar 2023
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 NOV 2020