



EXTENSION MEMO

Page 1 of 1

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


Catalog Number(s) / Size(s): K150ADJS-Series

Lot Number: K00S0142

Based on additional testing, the original date of expiration for this Kit has been extended by 12 months. The expiration date of K00S0142 has been extended to 31 AUG 2024.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
SARS-CoV-2 Spike Calibrator	A00S0046	31 AUG 2023	31 AUG 2024
Biotin SARS-CoV-2 Spike Antibody	B00T0064A	31 JUL 2024	31 AUG 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		15 SEP 2023

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



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K150ADJS-Series

Lot Number: K00S0142

Expiration Date: 31 AUG 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ SARS-CoV-2 Spike Antibody	D00T0064A	2-8°C	31 Aug 2024
Biotin SARS-CoV-2 Spike Antibody	B00T0064A	2-8°C	31 Jul 2024
SARS-CoV-2 Spike Calibrator	A00S0046	≤ -70°C	31 Aug 2023
Diluent 61	M0610004	≤ -10°C	31 Aug 2025
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 Aug 2025
Blocker S1	R0010824	≤ -10°C	30 Jun 2025
S-PLEX Enhance E1	Y0550012	≤ -10°C	31 Aug 2025
S-PLEX Enhance E2	Y0560013	≤ -10°C	31 Aug 2025
S-PLEX Enhance E3	Y0570013	≤ -70°C	31 Jul 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580017	≤ -70°C	31 Aug 2025
S-PLEX Detect D2	Y0590013	≤ -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:

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

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
SARS-CoV-2 Spike	175-2800	175	fg/mL	75% - 125%	< 25%	45.6	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
SARS-CoV-2 Spike	20X Calibrator	7520000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		12 MAY 2023