



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX SARS-CoV-2 Spike Kit
Lot Number: K00S0291
Expiration Date: 31 MAR 2026
Catalog Number: K150ADJS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin SARS-CoV-2 Spike Antibody	B00T0231	2-8°C	31 MAR 2028
TURBO-BOOST SARS-CoV-2 Spike Antibody	D00T0231	2-8°C	29 FEB 2028
Blocker S1	R0140008	≤ -10°C	31 JUL 2026
Diluent 61	M0610005	≤ -10°C	30 NOV 2027
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560017	≤ -10°C	31 JUL 2026
SARS-CoV-2 Spike Calibrator	A00S0116	≤ -70°C	31 MAR 2027
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580023	≤ -70°C	31 JUL 2026
S-PLEX Detect D2	Y0590019	≤ -70°C	30 SEP 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:

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

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
SARS-CoV-2 Spike	175 - 2800	700	fg/mL	75% - 125%	< 25%	28.9	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
SARS-CoV-2 Spike	20X Calibrator	7520000	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		09 SEP 2025